




High-throughput screening with nucleosome substrate identifies small-molecule inhibitors of the human histone lysine methyltransferase NSD2

Received for publication, June 3, 2018. Published, Papers in Press, June 26, 2018, DOI 10.1074/jbc.RA118.004274

Nathan P. Coussens[‡], Stephen C. Kales[‡], Mark J. Henderson[‡], Olivia W. Lee[‡], Kurumi Y. Horiuchi[§], Yuren Wang[§], Qing Chen[§], Ekaterina Kuznetsova[§], Jianghong Wu[§], Sirisha Chakka[‡], Dorian M. Cheff[‡], Ken Chih-Chien Cheng[‡], Paul Shinn[‡], Kyle R. Brimacombe[‡], Min Shen[‡], Anton Simeonov[‡], Madhu Lal-Nag[‡], Haiching Ma[§],  Ajit Jadhav[‡], and  Matthew D. Hall^{‡1}

From the [‡]National Center for Advancing Translational Sciences, National Institutes of Health, Rockville, Maryland 20850 and the [§]Reaction Biology Corporation, Malvern, Pennsylvania 19355

Edited by Joel Gottesfeld

The histone lysine methyltransferase nuclear receptor-binding SET domain protein 2 (NSD2, also known as WHSC1/MMSET) is an epigenetic modifier and is thought to play a driving role in oncogenesis. Both NSD2 overexpression and point mutations that increase its catalytic activity are associated with several human cancers. Although NSD2 is an attractive therapeutic target, no potent, selective, and bioactive small molecule inhibitors of NSD2 have been reported to date, possibly due to the challenges of developing high-throughput assays for NSD2. Here, to establish a platform for the discovery and development of selective NSD2 inhibitors, we optimized and implemented multiple assays. We performed quantitative high-throughput screening with full-length WT NSD2 and a nucleosome substrate against a diverse collection of bioactive small molecules comprising 16,251 compounds. We further interrogated 174 inhibitory compounds identified in the primary screen with orthogonal and counter assays and with activity assays based on the clinically relevant NSD2 variants E1099K and T1150A. We selected five confirmed inhibitors for follow-up, which included a radiolabeled validation assay, surface plasmon resonance studies, methyltransferase profiling, and histone methylation in cells. We found that all five NSD2 inhibitors bind the catalytic SET domain and one exhibited apparent activity in cells, validating the workflow and providing a template for identifying selective NSD2 inhibitors. In summary, we have established a robust discovery pipeline for identifying potent NSD2 inhibitors from small-molecule libraries.

Epigenetic modifiers are widely recognized as targets for therapeutic intervention, due to their critical roles in regulating gene expression and chromatin integrity in addition to their dysregulation in a range of human pathologies. In particular,

the nuclear receptor-binding SET domain (NSD)² family of histone lysine methyltransferase enzymes, NSD1, NSD2/WHSC1/MMSET, and NSD3/WHSC1L1, have all been implicated as cancer therapeutic targets (1). The catalytic SET domain (Suppressor of variegation, Enhancer of zeste, Trithorax) of NSD family enzymes catalyzes the mono- and di-methylation of the ϵ -amine of lysine 36 of histone H3 (H3K36), utilizing the methyl donor *S*-adenosyl-L-methionine (SAM) (Fig. 1A). The nonoverlapping roles of the NSD family of enzymes in normal physiology are attributed to a variety of additional domains, including a PWWP (proline-tryptophan-tryptophan-proline) domain, plant homeodomain, zinc finger domains, and a high-mobility group box domain (1).

NSD2 has been implicated as a therapeutic target for a variety of cancers. Because the gene is located within the Wolf-Hirschhorn syndrome critical region of chromosome 4, NSD2 is also known as Wolf-Hirschhorn syndrome candidate 1 (WHSC1) (2). NSD2 was first described as a gene dysregulated by the t(4;14)(p16.3;q32.3) translocation in ~15% of multiple myeloma (MM) cases and is called MMSET (2–4). The translocation results in a fusion transcript of NSD2 with the immunoglobulin heavy chain and increased NSD2 expression. The t(4;14) translocation can cause overexpression of both NSD2 and fibroblast growth factor receptor 3 (FGFR3) (2, 3). However, NSD2 is thought to be the primary oncogenic driver of the t(4;14)⁺ MM subtype because NSD2 is universally overexpressed, whereas FGFR3 is not expressed in ~30% of MM cases (4–6). The role of NSD2 in driving t(4;14)⁺ MM pathogenesis was supported by knockdown of NSD2 in MM t(4;14)⁺ cell lines, which led to reduced growth and tumorigenesis (7–11).

This work was supported by the National Center for Advancing Translational Sciences (NCATS) Division of Pre-Clinical Innovation Intramural Program. K. Y. H., Y. W., Q. C., E. K., J. W., and H. M. are employed by Reaction Biology. The content is solely the responsibility of the authors and does not necessarily represent the official views of the National Institutes of Health.

This article was selected as one of our Editors' Picks.

¹ To whom correspondence should be addressed. Tel.: 301-217-5727; E-mail: hallma@mail.nih.gov.

² The abbreviations used are: NSD, nuclear receptor-binding SET domain; qHTS, quantitative high-throughput screening; SAM, *S*-adenosyl-L-methionine; SPR, surface plasmon resonance; HTRF, homogeneous time-resolved fluorescence; DZNep, 3-deazaneplanocin A; SAH, *S*-adenosyl-L-homocysteine; MMSET, multiple myeloma SET domain; MM, multiple myeloma; CLL, chronic lymphocytic leukemia; CRC, concentration-response curve; TCEP, tris(2-carboxyethyl)phosphine; PMSF, phenylmethylsulfonyl fluoride; HBSS, Hanks' balanced salt solution; BisTris, 2-[bis(2-hydroxyethyl)amino]-2-(hydroxymethyl)propane-1,3-diol; Bicine, *N,N*-bis(2-hydroxyethyl)glycine; MTase-GloTM, Methyltransferase-Glo; S/B, signal-to-background; LPA, lysophosphatidic acid; NCATS, National Center for Advancing Translational Sciences; CV, coefficient of variation.

Conversely, transfection of $t(4;14)^{-}$ cells with NSD2 promotes tumorigenesis and oncogenic transformation of primary cells via elevated levels of dimethylated H3K36 (H3K36me2) (12). Numerous studies have linked increased expression of NSD2 with globally increased levels of H3K36me2 (9, 12–21). High expression of NSD2 protein has been demonstrated in many different human cancer types, including bladder, brain, gastrointestinal, lung, liver, ovary, skin, and uterus (18, 20, 22–28).

Notably, NSD2 is among the most frequently mutated genes in pediatric cancer genomes (29). The NSD2 SET domain variant, E1099K, was identified in both acute lymphoblastic leukemia tumors and cell lines with increased H3K36me2 that lack the $t(4;14)$ translocation (21, 30). Sequence results of >1,000 pediatric cancer genomes, representing 21 different cancers, revealed the E1099K variant in 14% of $t(12;21)$ ETV6-RUNX1 containing acute lymphoblastic leukemias (21). NSD2 is also among the most frequently mutated genes found in mantle cell lymphoma tumors, where both E1099K and T1150A variants are observed (31). The E1099K variant has also been reported in chronic lymphocytic leukemia (CLL) and lung and stomach cancers (32–35). Recombinant NSD2 E1099K showed higher *in vitro* activity compared with the WT enzyme (21). Ectopic expression of NSD2 E1099K induced H3K36me2 and promoted transformation, whereas knockdown of the enzyme reduced cell line proliferation and tumorigenesis (21).

Although NSD2 is an attractive therapeutic target, few small molecule inhibitors have been reported, and none demonstrate the desirable attributes of high-quality chemical probes (36). The compound LEM-06 (IC_{50} = 900 μ M) was discovered by virtual screening against an NSD2 homology model (37). The antiparasitic drug suramin inhibits NSD2 (IC_{50} = 0.3–21 μ M) but is a pan-inhibitor of methyltransferases (38, 39) as well as other enzymes (40). Likewise, the nonspecific histone lysine methyltransferase inhibitor chaetocin (IC_{50} = 3–6 μ M) showed similar inhibition of NSD1–3 (39). The natural product sinefungin is a close structural analog of SAM and a modest inhibitor of NSD2 (IC_{50} = 26–30 μ M) (41, 42). Structure–activity relationships have been reported for sinefungin analogs, the most potent of which inhibited the SET domains of NSD2 (IC_{50} = 1.8 μ M) and SETD2 (IC_{50} = 0.29 μ M) (41). Also, a peptide inhibitor of NSD2, PTD2 (IC_{50} = 3–22 μ M), has been reported that was derived from the histone H4 sequence (43).

A major challenge in screening for small molecule inhibitors is that native NSD2 requires nucleosomes as a substrate (17). Interestingly, the apparent specific activity of NSD2 is higher with HeLa-derived nucleosomes compared with recombinant nucleosomes, which has been attributed to unknown modifications of the native substrate (17). Thus, native nucleosomes purified from HeLa are likely a more physiologically relevant substrate than recombinant nucleosomes. Recombinant NSD2 does not act on peptides and is thus not amenable to the commonly adapted histone-derived peptide screening platforms. Here, we report biochemical assay development and pilot-scale screening using full-length recombinant WT, E1099K, and T1150A NSD2 enzymes with native HeLa nucleosomes as a substrate. Chemical libraries were screened in a three-dose-point quantitative high-throughput screening (qHTS) format (44) with the Methyltransferase-Glo (MTase-GloTM) biolumi-

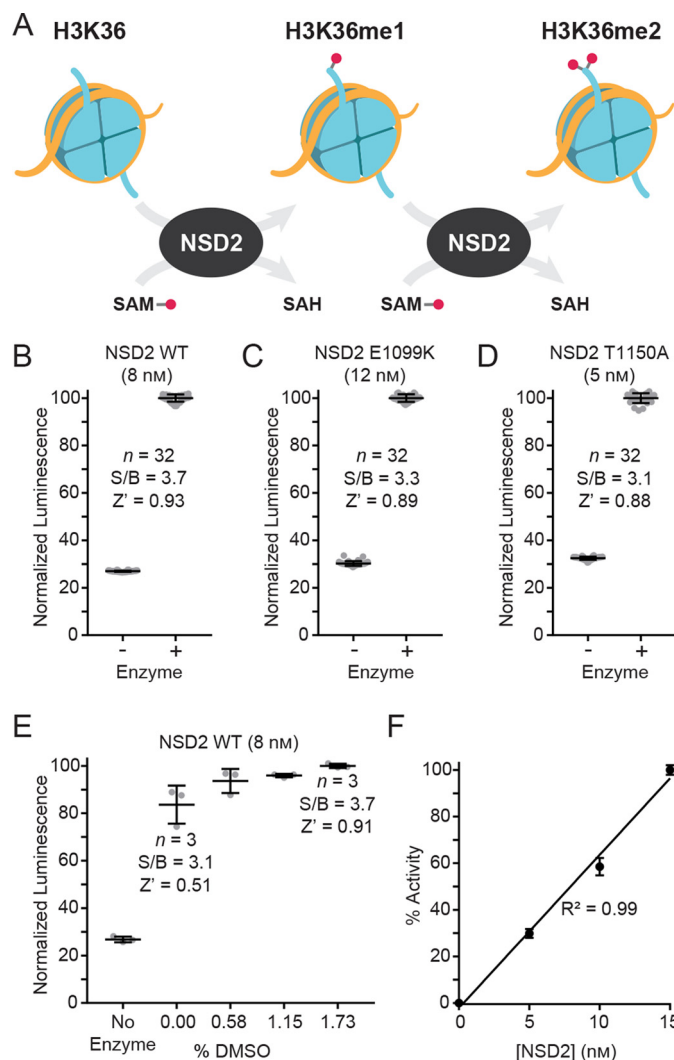


Figure 1. Optimization of the MTase-Glo primary assay. A, reaction scheme for the methylation of H3K36 by NSD2. B–D, optimized assay conditions for the full-length NSD2 WT (8 nM) (B), E1099K (12 nM) (C), or T1150A (5 nM) (D) enzymes with 500 nM nucleosomes, 0.58% DMSO, and 1 μ M SAM. The mean normalized luminescence values \pm S.D. ($n = 32$) are plotted for B–D. The assays are robust with S/B values >3 and Z' -factor values ~ 0.9 . E, titrations of DMSO demonstrated that the assay performance is not diminished by the introduction of vehicle up to 1.7% (mean \pm S.D.; $n = 3$). F, linear correlation is observed between the WT NSD2 enzyme concentration and methyltransferase activity (mean \pm S.D.; $n = 8$).

nescence assay (45). A number of counter and orthogonal assays were also developed to characterize the hits from the primary screen. Confirmed hits were validated with a radiolabeled SAM substrate-based NSD2 activity assay and further interrogated with binding studies by surface plasmon resonance (SPR), methyltransferase profiling, and activity assessments in U-2 OS cells.

Results

Primary assay development and optimization

To identify small molecule inhibitors of the WT NSD2, the MTase-Glo assay was optimized for use as a primary assay, with whole nucleosomes as the substrate, in 1,536-well format with 4- μ l reaction volumes (Fig. 1B). The MTase-Glo assay reagent measures methyltransferase activity through the coupling of

High-throughput NSD2 screening with nucleosome substrate

Table 1

Optimized protocol for the NSD2 MTase-Glo UltraGlo luciferase qHTS assay

Protocol Sequence for NSD2 MTase-Glo UltraGlo Luciferase qHTS Assay			
Step	Parameter	Value	Description
1	Reagent	4 μ l	SAH standard curve in reaction buffer [0 μ M - 1 μ M] as a 7-point 1:2 dilution series ($n = 4$), column 1
2	Reagent	3 μ l	Nucleosome [666.7 nM] in reaction buffer, columns 2-3
3	Reagent	3 μ l	NSD2 [10.7 nM] and nucleosome [666.7 nM] in reaction buffer, columns 4-48
4	Controls	23 nL	DMSO in columns 2-4
5	Library Compounds	23 nL	Columns 5-48
6	Time	30 min	Incubation
7	Reagent	1 μ l	S-adenosyl-L-methionine (SAM) [4 μ M], columns 2-48
8	Time	15 min	Incubation
9	Reagent	1 μ l	MTase-Glo Reagent [5X] in columns 1-48
10	Time	30 min	Incubation
11	Reagent	5 μ l	MTase-Glo Detection Solution in columns 1-48
12	Time	30 min	Incubation
13	Detection	Luminescence	ViewLux uHTS Microplate Imager (PerkinElmer)
Step	Notes		
1,2,3	White Medium Binding (MB) Greiner 1,536-well plates (789175-F); Reaction Buffer: 50 mM Tris-HCl, pH 8.8, 5 mM MgCl ₂ , 50 mM NaCl, 1 mM TCEP and 0.01% Tween 20		
4,5	Pintool transfer (tip wash sequence: DMSO, MeOH, 3 s vacuum dry)		
6, 8, 10 & 12	Room temperature		
7	Final reaction conditions (columns 4-48): 1 μ M SAM, 500 nM nucleosome, 8 nM NSD2, 50 mM Tris-HCl, pH 8.8, 5 mM MgCl ₂ , 50 mM NaCl, 1 mM TCEP and 0.01% Tween 20		
9	Conversion of SAH to ADP		
11	Conversion of ADP to ATP and detection by UltraGlo luciferase		
13	Settings: 30 s exposure, 1X binning		

the NSD2 reaction product S-adenosyl-L-homocysteine (SAH) to a bioluminescent signal (45). The assay was further optimized for the NSD2 variants E1099K (Fig. 1C) and T1150A (Fig. 1D), to enable insight into cross-inhibition of clinically-relevant NSD2 enzymes. The enzyme concentrations were optimized to allow a robust signal-to-background ratio (>3.0) in a 15-min reaction at room temperature while consuming $\leq 20\%$ substrate. The three optimized assays were robust with Z' -factor values near 0.9. There was no reduction in assay performance in the presence of up to 1.7% DMSO, which represents a triple aliquot of fixed volume (23 nL) transferred via a Kalypsys pintool equipped with a 1,536-pin array (Fig. 1E) (46). A titration of WT NSD2 demonstrated a linear correlation with methyltransferase activity as expected (Fig. 1F).

qHTS for NSD2 inhibitors

The primary screen was carried out by performing qHTS against 16,251 compounds at three concentrations (115, 57.5, and 11.5 μ M) in 1,536-well plates (Table 1). Commercially-available libraries screened included the LOPAC (1,280 compounds), Prestwick (1,360 compounds), MicroSource (2,000 compounds), and Tocris (1,304 compounds) collections. Addi-

tionally, four NCATS libraries were screened, including an epigenetics-focused collection (284 compounds), a natural products library (2,108 compounds), the NPACT library (5,099 compounds), and the NCATS Pharmaceutical Collection (2,816 compounds) (47). These libraries, enriched with pharmacologically active compounds, were selected to evaluate the suitability of the primary and secondary assays to identify NSD2 inhibitors. The overall quality of the primary screen data from 43 plates was high, with average values for the Z' -factor of 0.92 ± 0.02 , signal-to-background (S/B) of 3.28 ± 0.31 , and coefficient of variation (CV) for the DMSO control of $1.68 \pm 0.49\%$ (Fig. 2, A–C). The screen was conducted over a period of 4 months, so these data demonstrate excellent day-to-day reproducibility.

Hit confirmation and secondary screening

Of the 16,251 compounds evaluated in the primary screen, 536 compounds were active with a maximum response $\geq 50\%$, corresponding to a hit rate of 3.3% (Fig. 2D). Promiscuity scores were calculated for each hit, and any compounds with a promiscuity score higher than 0.2 were eliminated from the cherry-picking. Next, the remaining hits were evaluated by a structural

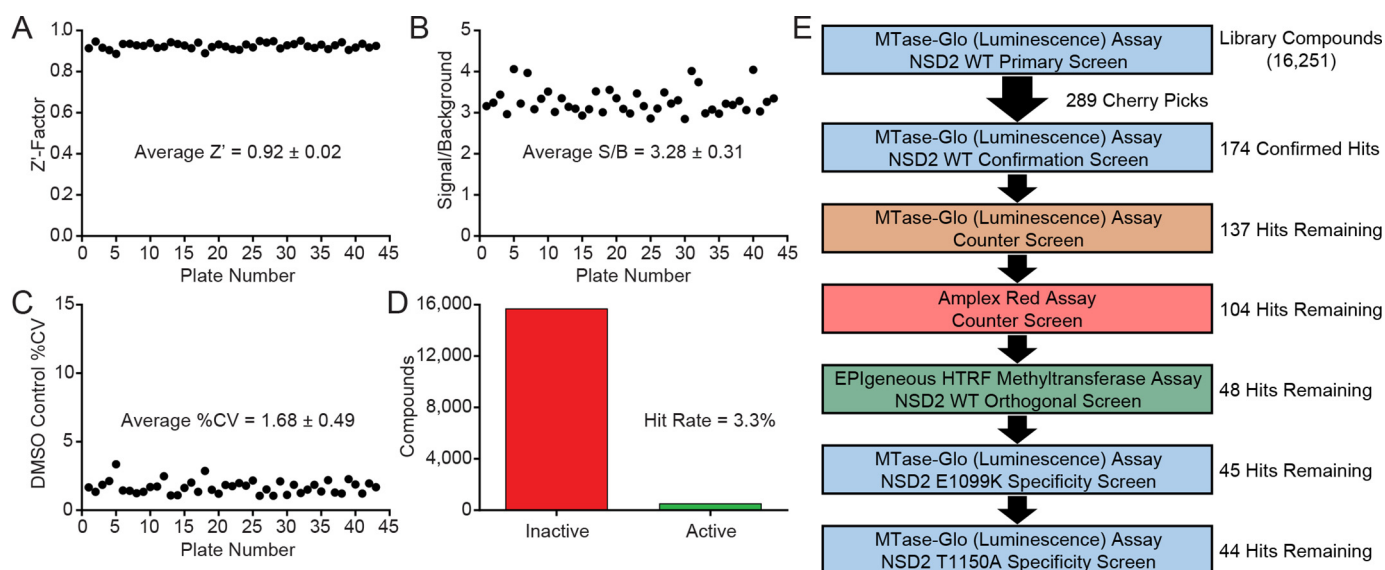


Figure 2. Quantitative high-throughput screening and secondary screening for inhibitors of NSD2 WT and variant enzymes. A–D, statistics from the qHTS primary screen against full-length WT NSD2 with the MTase-Glo methyltransferase assay for 43 1,536-well plates. A, overall assay performance of the primary screen was robust with an average Z'-factor value of 0.92. B, average signal-to-background ratio was 3.28. C, average % CV for the DMSO control was 1.68. D, of the 16,251 compounds screened, 536 were identified as active (3.3% hit rate). E, primary and secondary assay triage process. Eight compound libraries (16,251 compounds) were screened in qHTS format at three concentrations, resulting in 536 hits. From the initial hit list, 289 compounds were selected as cherry picks and further evaluated in 11-point dose response. Of the 289 cherry picks, the activities of 48 were confirmed against NSD2 with the primary and orthogonal assays with no activity observed by the counter assays. The majority of these compounds was also active against the NSD2 variants E1099K (45 compounds) and T1150A (44 compounds).

filter to eliminate the pan-assay interference compounds (PAINS) (48). Finally, 289 hit molecules were selected as cherry picks and prepared in 11-point concentration series from library stock solutions for further studies (Fig. 2E).

First, 174 of the cherry picks were confirmed as active in the primary assay (a 60% confirmation rate). Actives were defined as having concentration-response curves (CRCs) in the classes of 1.1, 1.2, 2.1, 2.2, and 3. In brief, classes 1.1 and 1.2 are the highest-confidence complete CRCs containing both upper and lower asymptotes with efficacies >80% and ≤80%, respectively. Classes 2.1 and 2.2 are incomplete CRCs having only one asymptote with efficacies >80% and ≤80%, respectively. Class 3 CRCs show activity at only the highest concentration or are poorly fit. Class 4 CRCs are inactive, having a curve-fit of insufficient efficacy or lacking a fit altogether (44). Second, an MTase-Glo counter screen was implemented without NSD2 but containing 200 nM SAH (mimicking 20% substrate conversion), which identified 37 assay interference compounds that might act by inhibiting the coupling enzymes, luciferase or the luminescent signal. Potential redox cycling by compounds was assessed with an Amplex Red assay performed in the presence of reducing agents (49), and 63 compounds were found to be active with a threshold of 3σ. To provide additional evidence for on-target activity against NSD2, the orthogonal EPIgenous homogeneous time-resolved fluorescence (HTRF) methyltransferase assay was utilized with reaction conditions identical to the primary assay (Fig. 3). The assay measures the NSD2 reaction product SAH, which competitively displaces d2-labeled SAH that is pre-bound to anti-SA-H labeled with Lumi4-Tb, resulting in a loss of FRET signal (50).

Among the 37 compounds found to be active with the MTase-Glo counter assay was a known luciferase inhibitor NCGC00183809 (Fig. 4A). The compound shares structural

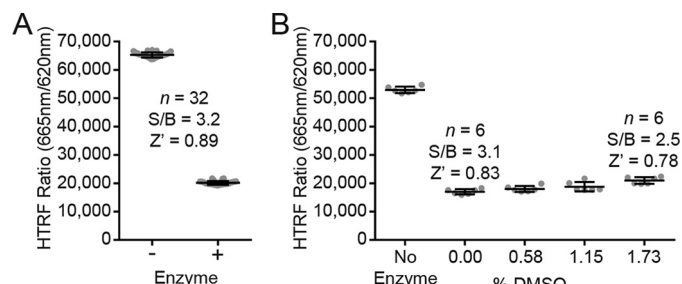


Figure 3. Optimized EPIgenous HTRF methyltransferase assay with the full-length NSD2 WT enzyme (8 nM), 500 nM nucleosomes, 0.58% DMSO, and 1 μM SAM. A, mean HTRF ratio values ± S.D. for n = 32 replicates are plotted, showing robust performance with S/B values >3 and a Z'-factor value of ~0.9. B, titrations of DMSO demonstrated that the assay performance is not diminished by the introduction of vehicle (mean ± S.D.; n = 6).

commonality with the firefly luciferase inhibitor PTC-124 (51, 52). As such, it was likely a false-positive hit that inhibits the UltraGlo luciferase enzyme utilized in the MTase-Glo assay. This notion was supported by the results of the orthogonal assay, which did not indicate inhibition of NSD2 activity. UltraGlo luciferase is a genetically evolved firefly luciferase containing 70 mutations to improve its robustness, thermal stability, and resistance to interference compounds (45, 53, 54), which is consistent with the 158-fold reduced potency of NCGC00183809 to UltraGlo compared with firefly luciferase (Fig. 4A). The nonspecific methyltransferase inhibitor 3-deazaneplanocin A (DZNep) showed similar activities in both the MTase-Glo primary and counter assays; however, no inhibition of NSD2 activity was observed with the orthogonal assay (Fig. 4B).

After the orthogonal assay, 48 confirmed hits remained that were not active in the two counter screens (Fig. 2E). Of the 48 confirmed inhibitors, 45 were also active against the E1099K

High-throughput NSD2 screening with nucleosome substrate

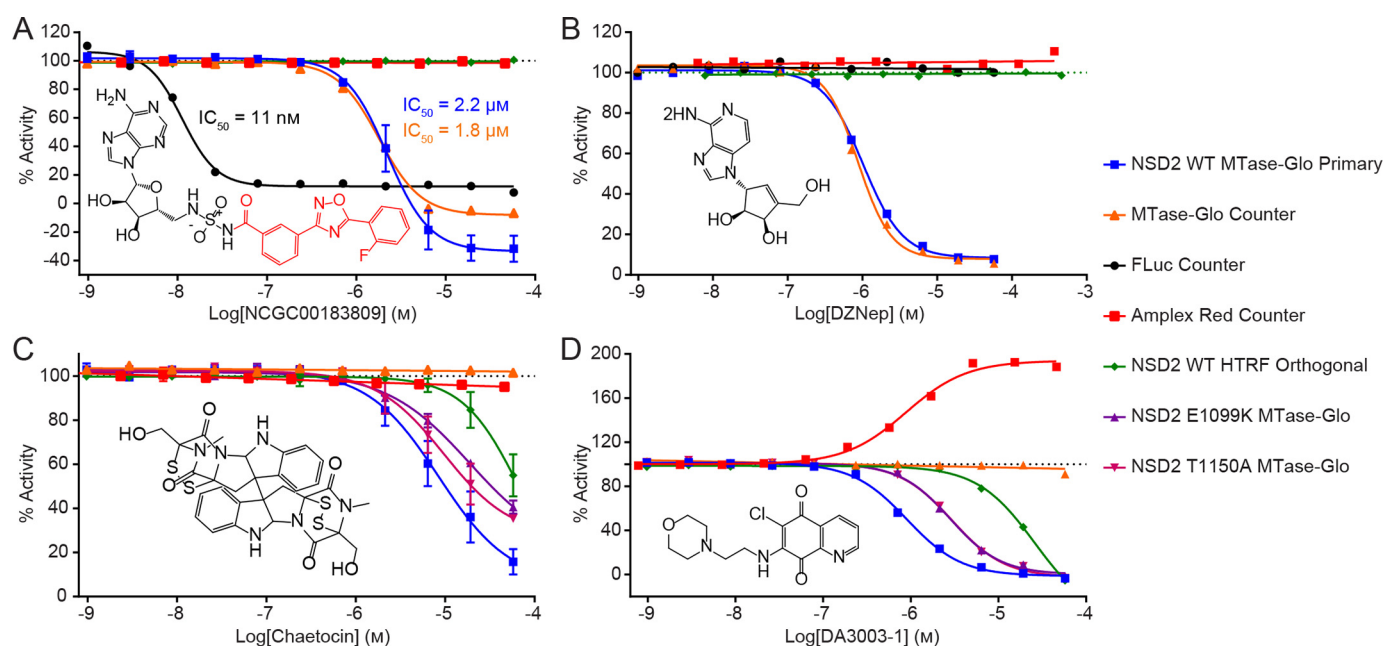


Figure 4. Primary and secondary assays characterize hit compounds against full-length NSD2 WT and variant enzymes. A, NCGC00183809 was identified by the primary assay ($IC_{50} = 2.2 \mu M$; Hill slope = -1.84) and contains the PTC-124 moiety (red) (51, 52). The compound showed similar activity with the counter assay ($IC_{50} = 1.8 \mu M$; Hill slope = -1.85) but was substantially more potent against firefly luciferase ($IC_{50} = 11 \text{ nM}$; Hill slope = -2.26). No activity against NSD2 was observed with the EPIgenous HTRF orthogonal assay. The data of WT NSD2 MTase-Glo are shown as the mean value \pm S.D. of $n = 2$ technical replicates. B, DZNep showed similar activities between the primary ($IC_{50} = 1 \mu M$; Hill slope = -1.63) and counter assays ($IC_{50} = 0.87 \mu M$; Hill slope = -1.95). No activity was observed with the orthogonal assay. C, confirmation of chaetocin as an NSD2 inhibitor. Chaetocin was identified as a hit from the primary screen, and the activity was confirmed by primary ($IC_{50} = 8.5 \mu M$; Hill slope = -1.09) and orthogonal assays ($IC_{50} = 67 \mu M$; Hill slope = -1.37). No activity was observed with the MTase-Glo counter assay or Amplex Red assay. Chaetocin also inhibits the NSD2 variants E1099K ($IC_{50} = 19 \mu M$; Hill slope = -0.93) and T1150A ($IC_{50} = 9.6 \mu M$; Hill slope = -1.15). Data from the MTase-Glo Counter and Amplex Red assays resulted from single experiments. All other data are shown as mean values \pm S.D. of n technical replicates ($n = 3$ for WT NSD2 MTase-Glo and $n = 2$ for HTRF orthogonal, NSD2 E1099K, and NSD2 T1150A). D, primary assay showed inhibition of WT ($IC_{50} = 0.9 \mu M$; Hill slope = -1.4), E1099K ($IC_{50} = 2.8 \mu M$; Hill slope = -1.51), and T1150A NSD2 ($IC_{50} = 2.9 \mu M$; Hill slope = -1.44) enzymes by DA3003-1, whereas no activity was observed with the counter assay. Inhibition of WT NSD2 was further supported by the orthogonal assay. However, DA3003-1 showed activity with the Amplex Red assay ($IC_{50} = 0.9 \mu M$; Hill slope = 1.25) indicating redox activity.

enzyme, and 44 showed activity against the T1150A variant. One of these hits is the nonspecific histone lysine methyltransferase inhibitor chaetocin, which was reported to inhibit NSD1, NSD2, and NSD3 (39). Chaetocin inhibited both WT and variant NSD2 enzymes without showing activity by the MTase-Glo counter assay or Amplex Red assay (Fig. 4C). The hit DA3003-1 is known to be redox-active (49, 55), and this was corroborated by the Amplex Red counter screen (Fig. 4D). Redox activity is undesirable, because it can result in nonspecific modulation of proteins, activation of cell pathways with redox-switches, and cytotoxicity (56). DA3003-1 was nevertheless selected for follow-up studies due to its submicromolar potency with the primary assay ($IC_{50} = 0.9 \mu M$). In addition to chaetocin and DA3003-1, three other hits were selected for further studies that have not been previously linked to NSD2 inhibition: PF-03882845, TC LPA5 4, and ABT-199 (Table 2). In anticipation of a full-fledged HTS, these five compounds were selected to validate our post-HTS workflow, which is intended to further evaluate compound activities.

Potency evaluation with the HotSpot radiolabel assay

Biochemical activities of the five selected hit compounds against WT and variant NSD2 enzymes were further validated by the radioisotope-based HotSpot assay (Fig. 5). The HotSpot assay incorporates [3H]SAM to assess total histone methylation by direct measurement of the filter-bound tritiated substrate, without the need for coupling enzymes or anti-

Table 2

Biochemical activities of five compounds that inhibit NSD2 WT and variant enzymes

All values are derived from one experiment unless followed by a symbol, in which case the value represents a global fit to n technical replicate experiments (*, $n = 2$; a , $n = 3$; b , $n = 4$; c , $n = 5$).

	Structure	DA3003-1	PF-03882845	Chaetocin	TC LPA5 4	ABT-199
Assay	Identifier	DA3003-1	PF-03882845	Chaetocin	TC LPA5 4	ABT-199
		NCGC00092289	NCGC00485023	NCGC00263643	NCGC00379183	NCGC00345789
MTase-Glo	NSD2	0.9	24	8.5 ^a	43	15
IC_{50} (μM)	WT					
	NSD2	2.8	34	19*	43	15
	E1099K					
	NSD2	2.9	38	9.6*	34	15
	T1150A					
HTRF	NSD2	26	31	67*	38	15
IC_{50} (μM)	WT					
HotSpot	NSD2	0.17*	7.6 ^a	0.13 ^a	8.5 ^a	1.7 ^a
IC_{50} (μM)	WT					
	NSD2	0.11	13	0.22	7.2	2.2
	E1099K					
	NSD2	0.17	11	0.6	8	2.3
	T1150A					
MTase-Glo Counter Assay		No Activity	No Activity	No Activity	No Activity	No Activity
IC_{50} (μM)						
Amplex Red Assay		0.9	No Activity	No Activity	No Activity	No Activity
IC_{50} (μM)						

bodies (38). Both DA3003-1 and chaetocin inhibited WT and variant NSD2 enzymes with submicromolar potencies, whereas PF-03882845, TC LPA5 4, and ABT-199 inhibited the enzymes at low-micromolar concentrations (Table 2).

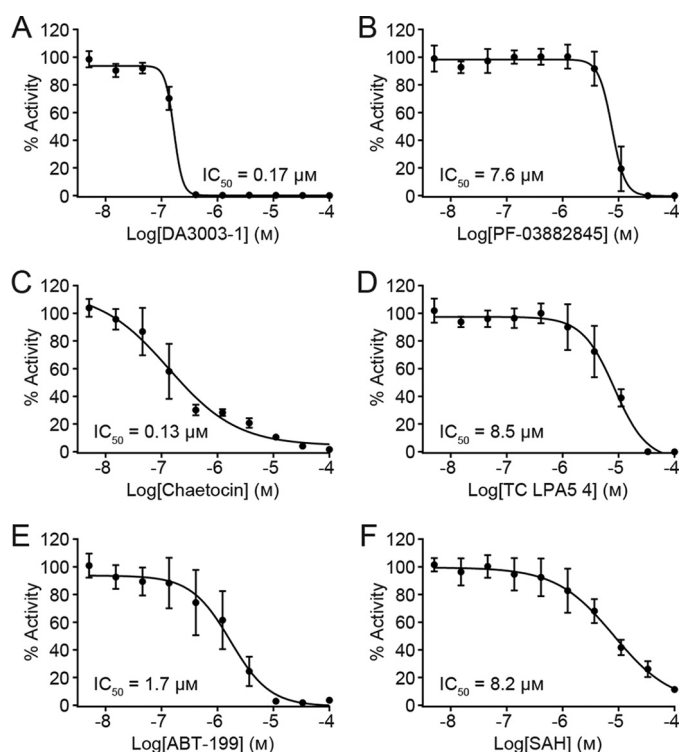


Figure 5. Inhibition of full-length WT NSD2 activity toward nucleosomes measured by the radiolabeled HotSpot assay with a 10-point concentration series of inhibitor. DA3003-1 potently inhibits NSD2 with an IC_{50} value of $0.17 \mu M$ (Hill slope = -5.57 ; $n = 4$) (A); PF-03882845 inhibits NSD2 activity with an IC_{50} value of $7.6 \mu M$ (Hill slope = -3.68 ; $n = 5$) (B); chaetocin is a potent inhibitor of NSD2 activity ($IC_{50} = 0.13 \mu M$; Hill slope = -0.71 ; $n = 5$) (C); TC LPA5 4 inhibits NSD2 with an IC_{50} value of $8.5 \mu M$ (Hill slope = -1.5 ; $n = 5$) (D); ABT-199 inhibits NSD2 with an IC_{50} value of $1.7 \mu M$ (Hill slope = -1.23 ; $n = 5$) (E). The positive control SAH inhibits NSD2 with an IC_{50} value of $8.2 \mu M$ (Hill slope = -0.85 ; $n = 4$) (F). Data are plotted as the mean value \pm S.D. of n technical replicate experiments.

Although chaetocin is known to inhibit NSD2, inhibition by DA3003-1, PF-03882845, TC LPA5 4, and ABT-199 has not been reported. Notably, the HotSpot assay was consistently more sensitive than the MTase-Glo and HTRF assays for both WT and variant enzymes (Table 2). Overall, the reaction conditions were very similar. The MTase-Glo and HTRF assays utilized 8 nM NSD2, 500 nM nucleosomes, 50 mM Tris-HCl, pH 8.8, 5 mM $MgCl_2$, 50 mM NaCl, 1 mM TCEP, 0.01% Tween; and the HotSpot assay used 10 nM NSD2, 400 nM nucleosomes, 50 mM Tris-HCl, pH 8.5, 5 mM $MgCl_2$, 50 mM NaCl, 1 mM DTT, 0.01% Brij35. A comparison of IC_{50} values determined by the HotSpot assay for all five compounds from reactions containing either 1 mM TCEP or DTT did not suggest that potency differences were due to the reducing agent (data not shown). Each of the five compounds inhibited WT, E1099K, and T1150A NSD2 enzymes with similar potencies (Table 2).

Direct binding of inhibitors to the NSD2 SET domain

To support an on-target mechanism of action for the five inhibitors, SPR was used to determine whether each inhibitor interacts with the catalytic SET domain of NSD2. As expected, the two positive controls, cofactor SAM and product SAH, both bound the NSD2 SET domain with low micromolar affinities (Table 3). With the exception of chaetocin, the inhibitors bound the NSD2 SET domain stoichiometrically, with dissoci-

Table 3

Direct binding of hit compounds to the NSD2 SET domain as measured by surface plasmon resonance, with SAH and SAM as positive controls

Compound name	k_{on} $M^{-1}s^{-1}$	k_{off} s^{-1}	K_d μM
SAH	2.33×10^3	1.19×10^{-2}	5.1
SAM	4.06×10^3	1.70×10^{-2}	4.2
DA03003-1	3.67×10^4	1.34×10^{-2}	0.37
PF-03882845	8.55×10^3	3.31×10^{-2}	3.9
Chaetocin	2.21×10^3	4.59×10^{-5}	0.02
TC LPA5 4	4.29×10^3	3.59×10^{-2}	8.4
ABT-199	4.39×10^3	3.64×10^{-2}	8.3

ation constants (K_d) comparable with the *in vitro* IC_{50} values determined by the HotSpot assay (Table 2). Chaetocin bound the SET domain with an apparent dissociation constant of 20 nM. Additionally, the data indicated super-stoichiometric binding, which might be due to a binding ratio higher than 1:1 that is consistent with chaetocin's two disulfide moieties forming adducts with the NSD2 protein or compound aggregation (data not shown) (57). These data indicate that all five compounds might mediate inhibition of NSD2 by directly binding to the catalytic SET domain.

Selectivity assessment by methyltransferase profiling

Methyltransferase profiling of the five NSD2 inhibitors was performed to examine the selectivity toward NSD2 compared with other methyltransferases. Activities of the inhibitors were evaluated against 35 other methyltransferases, including NSD1 and NSD3, with the HotSpot methyltransferase assay technology (Table 4). DA3003-1 was the least selective methyltransferase inhibitor and showed potent inhibition of nearly every enzyme except DOT1L (no activity) and GLP (weak inhibition). Furthermore, DA3003-1 inhibited 27 enzymes potently with submicromolar IC_{50} values. Both PF-03882845 and TC LPA5 4 inhibited most enzymes, albeit with weak potencies, but showed some selectivity to the PRMT5-MEP50 complex and the MLL4 complex. Chaetocin inhibited 18 methyltransferases, 12 with submicromolar potencies (including all NSD enzymes) and two with IC_{50} values above $75 \mu M$. ABT-199 inhibited 23 methyltransferases, two of which have IC_{50} values above $65 \mu M$. The NSDs were among the most potently inhibited enzymes, with IC_{50} values in the low-micromolar range.

Inhibitor activities in a cell-based assay

To evaluate the activities of the five NSD2 inhibitors in cells, U-2 OS human osteosarcoma cells were chosen due to a relatively high expression of endogenous NSD2 protein (18). To establish a baseline for reduced NSD2 activity, siRNA was used to knock down NSD2 in U-2 OS cells (Fig. 6A). Variations in NSD2 transcript knockdown were observed among the three unique siRNA molecules, and greater reductions in NSD2 corresponded to lower levels of H3K36me2 (Fig. 6, B and C). Knockdown of NSD2 did not appear to reduce the proliferation of U-2 OS cells (Fig. 6D). Next, the cells were treated with each of the five compounds in 10-point dose response from 0.0025 to $50 \mu M$ for 96 h. NSD2 inhibition should result in reduced H3K36me2. Total histone H3 and H3K36me2 levels were measured by Western blot analysis. The positive control DZNep was tested at a concentration of $10 \mu M$ in parallel with each

High-throughput NSD2 screening with nucleosome substrate

Table 4

Selectivity of inhibitor compounds toward WT NSD2 was assessed by profiling 37 other methyltransferases, including the E1099K and T1150A NSD2 variants, with the HotSpot methyltransferase assay

The IC₅₀ values for all test compounds are colored as a heat map with green indicating potent inhibition and red indicating weak inhibition. No color indicates no measurable activity. All values are derived from a single experiment except for those of WT NSD2, which result from a global fit of multiple data sets as indicated in Table 2.

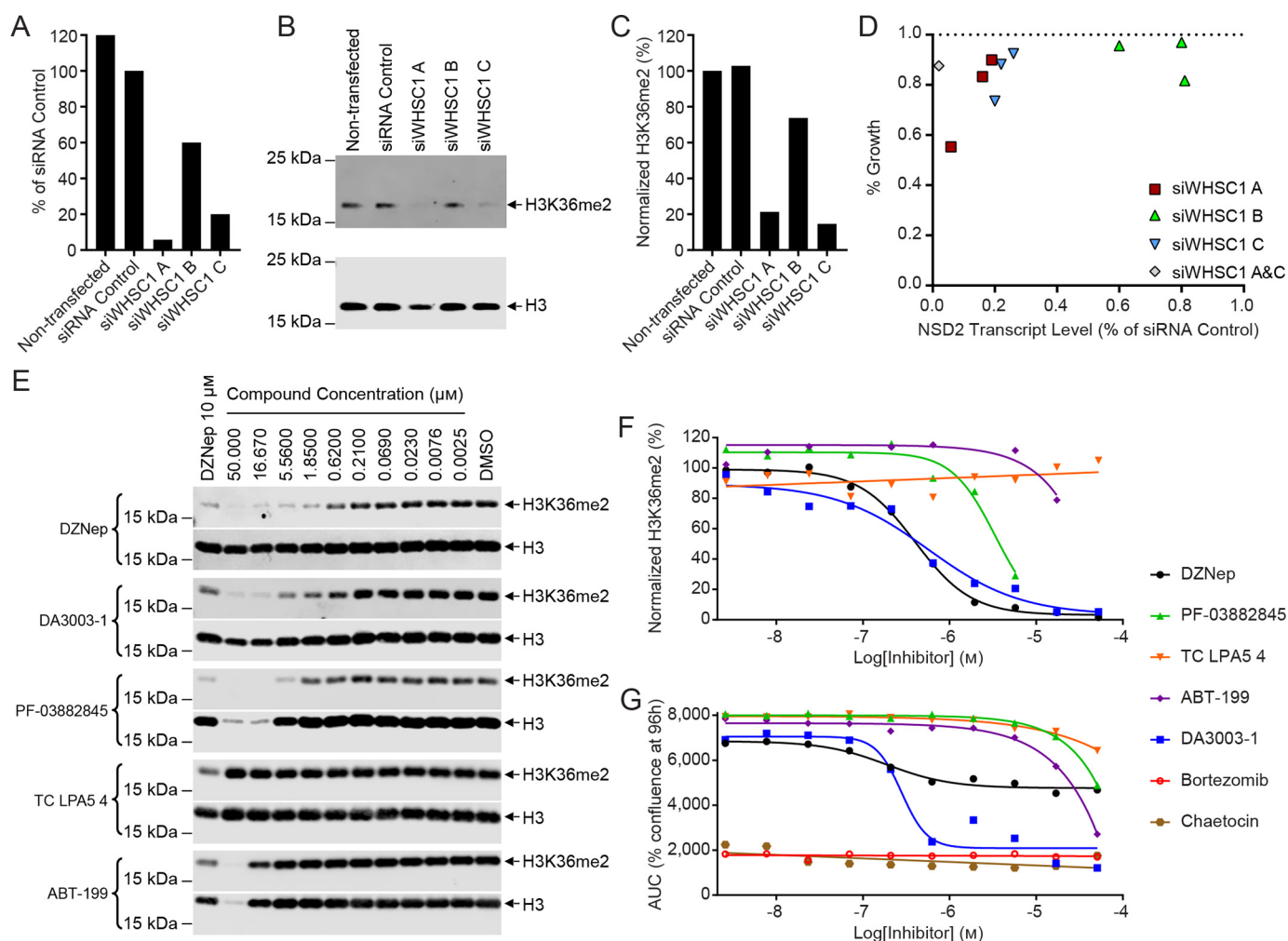
Methyltransferase	Substrate	DA3003-1 IC ₅₀ (μM)	PF-03882845 IC ₅₀ (μM)	Chaetocin IC ₅₀ (μM)	TC LPA5 4 IC ₅₀ (μM)	ABT-199 IC ₅₀ (μM)	Control IC ₅₀ (μM)	Control ID
ASH1L	Nucleosomes	0.2	9.8	0.31	9.2	3.3	0.1	Chaetocin
DNMT1	Poly (dl-dC)	1.7	14	No Inhibition	12	44	0.13	SAH
DNMT3a	Lambda DNA	1.1	18	No Inhibition	16	18	0.12	SAH
DNMT3b	Lambda DNA	0.38	19	No Inhibition	24	>100.00	0.065	SAH
DNMT3b-3L	Lambda DNA	0.72	31	No Inhibition	30	No Inhibition	0.012	SAH
DOT1L	Nucleosomes	No Inhibition	14	No Inhibition	8.2	2	0.74	SAH
EZH1 Complex	Core Histone	0.37	4	No Inhibition	40	66	19	SAH
EZH2 Complex	Core Histone	0.46	3.8	No Inhibition	40	23	17	SAH
EZH2 (Y641F) Complex	Core Histone	0.68	3.8	No Inhibition	31	21	68	SAH
G9a	Histone H3 (1-21)	0.9	No Inhibition	2.6	No Inhibition	No Inhibition	0.71	SAH
GLP	Histone H3 (1-21)	36	No Inhibition	77	No Inhibition	No Inhibition	0.53	SAH
MLL1 Complex	Nucleosomes	0.25	6.1	0.32	3.2	0.72	0.55	SAH
MLL2 Complex	Nucleosomes	0.32	15	3	18	4.6	27	SAH
MLL3 Complex	Core Histone	1.1	29	1.2	63	No Inhibition	57	SAH
MLL4 Complex	Nucleosomes	0.22	1.3	0.07	1	0.52	2.5	SAH
NSD1	Nucleosomes	0.11	12	0.06	12	2.4	9.7	SAH
NSD2	Nucleosomes	0.17	7.6	0.13	8.5	1.7	6.2	SAH
NSD2 (E1099K)	Nucleosomes	0.11	13	0.22	7.2	2.2	0.94	SAH
NSD2 (T1150A)	Nucleosomes	0.17	11	0.6	8	2.3	1.5	SAH
NSD3	Nucleosomes	0.26	3.2	0.33	2.3	1.3	0.09	Chaetocin
PRDM9	Histone H3	0.37	85	5.9	No Inhibition	No Inhibition	8.1	Chaetocin
PRMT1	Histone H4	0.26	8.5	No Inhibition	68	83	0.17	SAH
PRMT3	Histone H4	0.31	29	No Inhibition	5.7	36	2.3	SAH
PRMT4	Histone H3	1.4	29	No Inhibition	No Inhibition	No Inhibition	0.29	SAH
PRMT5	Histone H2A	0.65	11	No Inhibition	17	No Inhibition	49	SAH
PRMT5-MEP50 Complex	Histone H2A	0.21	0.38	No Inhibition	0.33	No Inhibition	1	SAH
PRMT6	GST-GAR	1	70	No Inhibition	No Inhibition	No Inhibition	0.084	SAH
PRMT7	GST-GAR	0.42	No Inhibition	156	No Inhibition	No Inhibition	0.045	SAH
PRMT8	Histone H4	0.22	21	No Inhibition	71	34	0.063	SAH
SET1b Complex	Core Histone	0.26	4.3	0.19	8	7.1	1.5	SAH
SET7/9	Core Histone	1.4	31	No Inhibition	52	No Inhibition	83	SAH
SET8	Nucleosomes	4.4	42	No Inhibition	44	5.3	0.95	Ryuvidine
SETD2	Nucleosomes	0.21	13	0.45	13	3.3	5.2	SAH
SMYD1	Core Histone	1	11	No Inhibition	21	No Inhibition	37	SAH
SMYD2	Histone H4	1.7	32	No Inhibition	21	44	4.5	LLY 507
SUV39H1	Histone H3	0.47	No Inhibition	0.74	No Inhibition	No Inhibition	25	SAH
SUV39H2	Histone H3	0.62	No Inhibition	0.57	No Inhibition	No Inhibition	55	SAH
SUV420H1TV2	Nucleosomes	0.34	4.4	No Inhibition	4.2	2.2	31	SAH

compound and was also evaluated in dose response (Fig. 6E). Densitometry was used to quantify both H3K36me2 and total H3, and the densities of H3K36me2 were normalized to those of H3 (Fig. 6F). The growth of U-2 OS cells over 96 h in the presence of test compounds at the same concentrations was also evaluated (Fig. 6G). The IC₅₀ value of DZNep for reducing H3K36me2 levels in U-2 OS cells was 390 nM with a modest dose-dependent reduction in U-2 OS confluency (IC₅₀ = 180 nM). Similar to DZNep, DA3003-1 treatment also resulted in a dose-dependent reduction in H3K36me2 (IC₅₀ = 545 nM); however, higher drug concentrations resulted in cytotoxicity (CC₅₀ = 270 nM) similar to chaetocin and the control bortezomib, both of which were cytotoxic at all concentrations tested. PF-03882845 reduced H3K36me2 (IC₅₀ = 3.2 μM) over

a range of concentrations that had negligible influence on growth; however, substantial cell death at concentrations above 16.7 μM was observed during the Western blot analysis experiment (Fig. 6E), so data from the two highest concentrations were excluded from the potency calculation. TC LPA5 4 did not reduce H3K36me2 over the concentrations tested, and minimal reduction in growth was observed. ABT-199 modestly reduced H3K36me2 levels at 16.7 μM; however, higher concentrations were cytotoxic.

Discussion

A number of histone lysine methyltransferases have been implicated as attractive therapeutic targets in the field of oncology, and small molecule inhibitors are in different stages of



preclinical and clinical development (59, 60). Selective inhibitors of NSD2 are of major interest both to advance basic research and for therapeutic development. However, NSD2 is regarded as a challenging target (61), and no selective small molecule inhibitors of NSD2 have been reported to date. Challenges in studying NSD2 *in vitro* include that the target of NSD2 methylation depends on the nature of the substrate (17), and full-length NSD2 is only active against a nucleosome substrate (38), which may be cost prohibitive in many cases.

A wide variety of methyltransferase assays have been described in the literature (38, 45, 62–67); however, few optimized assays have been reported for NSD2 in plate formats to enable HTS. Recently, a radiolabeled [3 H]SAM assay was optimized for the NSD2 SET domain with both histone octamer (Z' = 0.69) and nucleosome (Z' = 0.75) substrates in a 384-well format (39). The assay utilizing a nucleosome substrate was

used to screen 1,040 compounds from the Prestwick Chemical Library at a single concentration of 25 μ M. Although the reported assays show robust performance in the 384-well format, the use of radiolabeled reagents for HTS is a challenge for many laboratories due to safety regulations and disposal costs.

Herein, we report the implementation and validation of two optimized homogeneous NSD2 activity assays in the highly miniaturized 1,536-well format for the identification of small molecule inhibitors from chemical libraries. The assays were utilized to screen the full-length WT NSD2 enzyme against a nucleosome substrate in qHTS format with three concentrations of test compounds. The use of qHTS reduces both false-positive and false-negative hits common to single-point HTS and facilitates selection of actives (44). For the primary screen, we used the recently reported Methyltransferase-Glo assay reagent with a sensitive biolumines-

cent readout (45). A similar approach has been applied for the discovery of NSD1 inhibitors by HTS (68). After screening eight libraries, including numerous pharmacologically active collections, containing over 16,000 compounds, many hits were identified, including chaetocin, which is known to inhibit NSD2 (39). By incorporating orthogonal and counter screens, hits were prioritized for subsequent follow-up studies.

Among the confirmed active inhibitors, DA3003-1, PF-03882845, chaetocin, TC LPA5 4, and ABT-199 were selected for further characterization. *In vitro* potencies were determined by the HotSpot assay, which is a direct readout of the NSD2 reaction product (38). The HotSpot assay is very similar in format to the traditional gold standard radioisotope detection used in conjunction with gel electrophoresis or MS. The five compounds inhibited WT NSD2 as well as the E1099K and T1150A variant enzymes.

DA3003-1 is a cell-permeable Cdc25 phosphatase inhibitor that potently and irreversibly inhibits all Cdc25 isoforms, including Cdc25A ($IC_{50} = 29$ nM), Cdc25B2 ($IC_{50} = 95$ nM), and Cdc25C ($IC_{50} = 89$ nM) (69). It is known that DA3003-1 is capable of redox cycling (49, 55), which was verified here by the Amplex Red counter screen. In addition to potent inhibition of NSD2 activity, our data demonstrate that DA3003-1 bound the SET domain with a strong affinity ($K_d = 370$ nM, Table 3), which is comparable with its potency ($IC_{50} = 170$ nM, Table 2). Together, this suggests that DA3003-1 inhibits NSD2 through a direct interaction with the catalytic SET domain, although it is most likely nonspecific. Notably, DA3003-1 inhibited 27 of the 38 methyltransferase enzymes tested with submicromolar potencies and another eight with IC_{50} values near 1 μ M. The potency of DA3003-1 in cells with respect to reducing H3K36me2 ($IC_{50} = 545$ nM) corresponds to cytotoxicity ($CC_{50} = 270$ nM).

The Pfizer compound PF-03882845 is a highly potent mineralocorticoid receptor antagonist ($IC_{50} = 0.755$ nM) (70). In comparison, it inhibited NSD2 *in vitro* with an $\sim 10,000$ -fold weaker potency ($IC_{50} = 7.6$ μ M) and bound the SET domain ($K_d = 3.9$ μ M) within 2-fold of the IC_{50} value. Profiling studies indicated that PF-03882845 inhibits many other methyltransferases with modest potencies, although it inhibits the histone arginine methyltransferase PRMT5 with a submicromolar potency. Interestingly, PF-03882845 reduced H3K36me2 in cells with a potency (3.2 μ M) within 2-fold of the biochemical IC_{50} value. Concentrations at and above 17 μ M resulted in substantial cell death in the Western blot analysis experiment, but only minimal reductions in cell growth were observed with up to 50 μ M drug at 96 h in the cell growth assay.

TC LPA5 4 was first reported by Sanofi Aventis as a selective lysophosphatidic acid receptor 5 (LPA₅) antagonist that inhibited LPA-mediated human platelet aggregation with an IC_{50} value of 2.2 μ M (71). The biochemical potency of TC LPA5 4 against NSD2 ($IC_{50} = 8.5$ μ M) was nearly identical to its affinity to the catalytic SET domain ($K_d = 8.4$ μ M). The methyltransferase activity profiles of TC LPA5 4 and PF-03882845 showed striking similarities (Table 4). Both compounds inhibited WT NSD2 with an IC_{50} value near 8 μ M. Also, both compounds were most potent against the PRMT5–MEP50 complex with

nearly identical IC_{50} values. Furthermore, both compounds had ~ 3 -fold weaker potency against the MLL4 complex. Compared with the PRMT5–MEP50 complex, the potency values of PF-03882845 were at least 8-fold weaker against 31 other methyltransferases. Similarly, compared with the PRMT5–MEP50 complex, the potency value of TC LPA5 4 is at least 7-fold weaker against 28 other methyltransferases. No activity or toxicity was observed for TC LPA5 4 in U-2 OS cells.

The fungal mycotoxin chaetocin is known to inhibit NSD2 (39), so identifying it as a hit further validated our screening approach. Of the 38 methyltransferases profiled, chaetocin inhibited 12 with submicromolar potencies. Notably, the methyltransferase profiling indicated chaetocin potencies of 740 nM against SUV39H1 and 570 nM against SUV39H2, which is consistent with a previously reported value of 600 nM (72). Chaetocin was initially reported to be a specific inhibitor of the histone lysine methyltransferase SU(VAR)3-9 both *in vitro* and *in vivo* (72). However, the two disulfide bonds of chaetocin can complicate bioassay interpretation because of the potential for redox activity and covalent modification of proteins (36). Indeed, reports have indicated that the activity against histone lysine methyltransferases is due to chemical modification of the enzyme by the disulfide groups (73, 74). Superstoichiometric binding of chaetocin to the NSD2 SET domain was observed by surface plasmon resonance, which might be due to the formation of direct compound–thiol adducts. The affinity of chaetocin to the NSD2 SET domain was strong ($K_d = 20$ nM) with a biochemical potency about 7-fold weaker. Chaetocin was cytotoxic to U-2 OS cells at all concentrations tested.

ABT-199, also known as GDC-0199 or venetoclax, binds BCL-2 with a subnanomolar affinity ($K_i < 0.010$ nM) and is approved by the Food and Drug Administration for the treatment of CLL (75, 76). The biochemical potency of ABT-199 against WT NSD2 ($IC_{50} = 1.7$ μ M) was similar against the NSD2 variants, NSD1 and NSD3 (Table 4). The compound was most potent against the MLL1 and MLL4 complexes. The affinity of ABT-199 to the NSD2 SET domain ($K_d = 8.3$ μ M) is nearly 5-fold weaker than the biochemical potency. At a concentration of 50 μ M, ABT-199 was cytotoxic, consistent with its use in oncology, but any influences on cellular H3K36me2 levels were negligible.

DZNep is a carbocyclic analog of adenosine and a derivative of the antibiotic neplanocin-A. It was initially reported as a competitive inhibitor of SAH hydrolase at picomolar concentrations (77). DZNep has since been reported as a global histone methylation inhibitor when used at substantially higher concentrations (36, 78, 79). Although DZNep was identified as a hit from the primary screen, similar activities were observed with the MTase-Glo primary and counter assays, and no inhibition of NSD2 activity was observed with the orthogonal assay. The potency of DZNep in reducing H3K36me2 in U-2 OS cells ($IC_{50} = 390$ nM; Fig. 6F) is similar to its potency in reducing H3K27me3 in SU-DHL-6 cells ($IC_{50} = 160$ nM) when assessed by the same method of Western blot analysis (80). DZNep has been also shown to reduce H3K36me2 in SW480 cells at a concentration of 5 μ M (81).

The purpose of this study was to establish an HTS discovery pipeline for NSD2 and to evaluate the workflow for identifying

high-quality tool inhibitors of NSD2. The majority of the molecules screened were from pharmacologically active libraries that served to validate the primary and secondary assays. The identification of known methyltransferase inhibitors, including chaetocin and DZNep, further validated the workflow. During this pilot, many known interference compounds were identified by the secondary assays, thereby demonstrating how such bad actors behave among the various assays. Five actives selected from the primary screen were shown to bind the catalytic SET domain and inhibit NSD2 activity *in vitro*. Although these studies confirm inhibition of NSD2, they do not rule out inhibition by intractable mechanisms of action, such as nonspecific reactivity, redox, or aggregation. Two of the five compounds reduced H3K36me2 in U-2 OS cells, but the mechanisms are likely to be complicated and involve multiple targets. These studies provide a basis for the future discovery and development of novel selective NSD2 inhibitors by establishing a robust workflow for identifying and triaging hits from high-throughput screens.

Experimental procedures

Chemicals, reagents, and libraries

MTase-Glo (V7602) was purchased from Promega, Inc. (Madison, WI). The EPIgeneous HTRF methyltransferase assay (catalog no. 62SAHPEB) was purchased from Cisbio (Bedford, MA). Dimethyl sulfoxide (DMSO) was purchased from ThermoFisher Scientific (Pittsburgh, PA). The libraries screened include the following: Library of 1280 Pharmacologically Active Compounds (LOPAC¹²⁸⁰; Sigma); Tocris (1,304 compounds, Tocris Bioscience, Bristol, UK); Prestwick (1,360 compounds, Prestwick Chemical, San Diego); MicroSource (2,000 compounds, MicroSource Discovery Systems, Gaylordville, CT); NPC (the NCATS Pharmaceutical Collection (2,816 compounds) (47); NPACT (NCATS Pharmacologically Active Chemical Toolbox, 5,099 compounds); an epigenetic collection (284 compounds); and a natural products library (2,108 compounds). The reference compound DZNep was purchased from Selleckchem and dissolved with DMSO to a 10 mM stock. The U-2 OS human osteosarcoma cell line and McCoy's 5A medium were purchased from American Type Culture Collection (Manassas, VA). U-2 OS cells were grown in McCoy's 5A medium supplemented with 10% fetal bovine serum (FBS), 100 µg/ml penicillin, and 100 µg/ml streptomycin. Cultures were maintained at 37 °C in a humidified atmosphere of 5% CO₂ and 95% air. The 12% BisTris gel, nitrocellulose membrane, Lipofectamine RNAiMAX reagent, and High-Capacity RNA-to-cDNA kit were purchased from ThermoFisher Scientific. The RNeasy mini kit and AllStars Hs Cell Death Control siRNA were obtained from Qiagen (Germantown, MD). The TaqMan Gene Expression Master Mix and TaqMan WHSC1 human (Hs00983720_m1) were obtained from Applied Biosystems (Foster City, CA). Silencer Select Negative Control No. 2 siRNA, siWHSC1 A (Assay ID 42800), siWHSC1 B (Assay ID 42732), siWHSC1 C (Assay ID 42650), and human GAPDH endogenous control (4352934E) were obtained from Life Technologies, Inc. Dimethyl histone H3 (Lys-36) rabbit mAb (2901) and histone H3 mouse mAb (3638) were purchased from Cell

Signaling Technology. Anti-rabbit IgG IRDye 680RD and anti-mouse IgG IRDye 800CW secondary antibodies were purchased from LI-COR.

Enzymes and substrates

Human ASH1L (residues 2046–2330; GenBankTM accession no. NM_018489) was expressed in *Escherichia coli* with an N-terminal polyhistidine tag. Human DNMT1 (residues 2–1632; NM_001130823) was expressed in an insect cell/baculovirus expression system as an N-terminal GST fusion. Human DNMT3a (residues 2–912; NM_175629) was expressed in an insect cell/baculovirus expression system as an N-terminal GST fusion. Human DNMT3b (residues 2–853; NM_006892) was expressed in an insect cell/baculovirus expression system as an N-terminal GST fusion. The two proteins human DNMT3b (residues 564–853; NM_006892) containing an N-terminal polyhistidine tag and human DNMT3L (residues 160–387; NM_013369) with an N-terminal GST tag were coexpressed in an insect cell/baculovirus expression system. Human DOT1L (residues 1–416; NM_032482) was expressed in *E. coli* as an N-terminal GST fusion. Human recombinant EZH1 (residues 2–747; NM_001991) or EZH2 (residues 2–746; NM_001203247) was coexpressed with human recombinants AEBP2 (2–517; NM_001114176), EED (2–441; NM_003797), RbAp48 (2–425; NM_005610), and SUZ12 (2–739; NM_015355) in an insect cell/baculovirus expression system to form the five-member EZH1 or EZH2 complexes. All proteins were full-length (residue 2 through C terminus). The EED subunit incorporated an N-terminal FLAG tag, and all others included an N-terminal polyhistidine tag. Human GLP (residues 894–1298; NM_024757) and human G9a (residues 786–1210; NM_006709) were expressed as N-terminal GST fusion proteins in *E. coli*. Human MLL1 (residues 3745–3969; NM_005933), human WDR5 (22–334; NM_017588), RbBP5 (1–538; NM_005057), Ash2L (2–534; NM_001105214), and DPY-30 (1–99; NM_0325742) were expressed in *E. coli* with N-terminal polyhistidine tags, assembled as a complex (two copies of DPY-30), and stored in 20 mM Tris-HCl, pH 7.5, 300 mM NaCl, 1 mM TCEP, 10% (w/v) glycerol, and 1 mM ZnCl₂. Human MLL2 (residues 5319–5537; NM_003482), human MLL3 (residues 4689–4911; NM_170606), and human MLL4 (residues 2490–2715; NM_014727) were expressed in *E. coli* with N-terminal polyhistidine tags, and SET1B (residues 1629–1923; NM_015048) was expressed in *E. coli* with an N-terminal GST tag. All four were assembled in complexes as MLL1, as described above. Human recombinant NSD1 (residues 1538–2696; NM_022455) was expressed with an N-terminal polyhistidine tag in an insect cell/baculovirus expression system. Human recombinant NSD2 (residues 2–1365; NM_001042424) was expressed with an N-terminal polyhistidine tag in an insect cell/baculovirus expression system. Human recombinant NSD2-SET domain for SPR studies (residues 934–1241; NM_001042424) was expressed with an N-terminal polyhistidine tag in *E. coli*. Human recombinant NSD3 (residues 1021–1322; NM_023034) was expressed in *E. coli* as an N-terminal GST fusion. Human PRDM9 (residues 2–414; NM_020227) was expressed in an insect cell/baculovirus expression system as an N-terminal GST fusion. Human

recombinant PRMT1 (residues 2–371; NM_001536) was expressed in *E. coli* as an N-terminal GST fusion. Human recombinant PRMT3 (residues 2–531; NM_005788) was expressed in *E. coli* with an N-terminal polyhistidine tag. Human recombinant PRMT4 (residues 2–608; NM_199141) was expressed in *E. coli* as an N-terminal GST fusion. Human recombinant PRMT5 (residues 2–637; NM_006109) was expressed with an N-terminal FLAG tag in an insect cell/baculovirus expression system. The complex of PRMT5–MEP50 was coexpressed with PRMT5 as described above with MEP50 (residues 2–342; NM_024102) containing an N-terminal polyhistidine tag in an insect cell/baculovirus expression system. Human recombinant PRMT6 (residues 2–375; NM_018137) and PRMT7 (residues 2–692; NM_019023.2) were expressed with N-terminal polyhistidine tags in an insect cell/baculovirus expression system. Human recombinant PRMT8 (residues 61–394; NM_019854) was expressed in *E. coli* with N- and C-terminal polyhistidine tags. Human recombinant SET7/9 (residues 2–366; NM_030648) was expressed in *E. coli* as an N-terminal GST fusion with a C-terminal polyhistidine tag. Human recombinant SETD2 (residues 1434–1711; NM_014159) and SMYD1 (residues 2–490; NM_198274), both with N-terminal GST fusions, were expressed in *E. coli*. Human recombinant SMYD2 (residues 2–433; NM_020197) was expressed in *E. coli* with an N-terminal polyhistidine tag. Human recombinant SUV39H1 (residues 44–412; NM_003173) and SUV39H2 (residues 48–410; NM_001193424), both with C-terminal polyhistidine tags, were expressed in *E. coli*. Human recombinant SUV420H1-tv2 (transcript variant 2, residues 2–393; NM_016028) was expressed as an N-terminal GST fusion in an insect cell/baculovirus expression system. Purified nucleosomes (HMT(35–130)) were obtained from HeLa according to Schnitzler (82). Core histones, including histone 5 (HMT(35–435)), were purified from chicken erythrocytes by a modification of the method of Schnitzler (82) followed by acid extraction/dialysis (83). Human recombinant histone H2A (residues 1–130; NM_021052) and histone H3.3 (residues 1–136; NM_002107), both untagged, were expressed in *E. coli*. Human recombinant GST-GAR (glycine- and arginine-rich sequence from the N terminus of fibrillarin, residues 2–78; NM_001436) was expressed in *E. coli* as an N-terminal GST fusion. The following substrates were purchased from vendors: poly(dI-dC)(dI-dC) from Sigma; λ DNA from New England Biolabs; histone H4 from BPS Biosciences, and histone H3(1–21) peptide from AnaSpec. The following reference compounds were purchased from vendors: SAH and chaetocin from Cayman Chemicals, and LLY 507 and ryludine from R & D Systems.

Methyltransferase-Glo assay

MTase-Glo assays were performed by a multistep format in white solid bottom 1,536-well plates (Greiner, catalog no. 789175-F). First, 23 nl of compounds (or DMSO control) were pin-transferred into 3 μ l of reaction buffer (50 mM Tris-HCl, pH 8.8, 5 mM MgCl₂, 50 mM NaCl, 1 mM TCEP, and 0.01% Tween 20) containing 666.7 nM (500 nM final) nucleosomes and either 10.7 nM (8 nM final) WT, 16 nM (12 nM final) E1099K, or 6.67 nM (5 nM final) T1150A NSD2 enzyme or no enzyme (low

activity control, columns 2–3). Plates were then incubated for 30 min at room temperature prior to reaction initiation with 1 μ l of 4 μ M (1 μ M final) SAM in reaction buffer and incubated at room temperature for 15 min. Upon completion, methyltransferase conversion of SAM to SAH was then detected using a two-step detection system where 1 μ l of MTase-Glo reagent was first added to each well to convert SAH to ADP for 30 min at room temperature. Secondly, 5 μ l of MTase-Glo Detection Solution was added to each well and allowed to incubate for 30 min at room temperature to convert ADP to ATP, which was then measured by luminescence detection using a ViewLux uHTS Microplate Imager (PerkinElmer Life Sciences) and compared with control samples to determine relative activity.

Methyltransferase-Glo counter assay

MTase-Glo counter assay was performed with identical procedures as with the MTase-Glo primary assay but without NSD2 enzyme or nucleosomes. Instead, 200 nM SAH was added to mimic the reaction with 20% substrate conversion.

EPiGenetic HTRF methyltransferase assay

The EPiGenetic HTRF methyltransferase assay was performed using white solid bottom 1,536-well plates (Greiner, catalog no. 789175-F). First, 23 nl of compounds (or DMSO control) were pin-transferred into 3 μ l of reaction buffer (50 mM Tris-HCl, pH 8.8, 5 mM MgCl₂, 50 mM NaCl, 1 mM TCEP and 0.01% Tween 20) containing 666.7 nM (500 nM final) nucleosomes and either 10.7 nM (8 nM final) WT NSD2 enzyme or no enzyme (low activity control, columns 2–3). Plates were then incubated for 30 min at room temperature prior to reaction initiation with 1 μ l of 4 μ M (1 μ M final) SAM in reaction buffer and incubated at room temperature for 15 min. After the incubation period, 0.8 μ l of EPiGenetic Detection Buffer One was added to each well, followed by a 10-min incubation at room temperature. Next, anti-SA-H/Lumi4-Tb solution was prepared according to the manufacturer's instructions, and 1.6 μ l of the solution were added to each well. Finally, the SAH-d2 conjugate was prepared as a 32-fold dilution according to the manufacturer's instructions, and 1.6 μ l of the solution was added to each well. The assay plate was allowed to incubate for 1 h at room temperature before detection of the HTRF signal using an Envision plate reader (PerkinElmer Life Sciences).

Amplex red (10-acetyl-3,7-dihydroxyphenoxazine) assay

The assay was adapted from a previously described protocol to assess redox cycling of compounds in the presence of reducing agents (49). 23 nl of compounds (or DMSO control) were pin-transferred into 2.5 μ l of HBSS (ThermoFisher Scientific; containing 1.26 mM CaCl₂, 0.49 mM MgCl₂, 1 g/liter D-glucose) in black 1,536-well plates. Compound fluorescence was measured immediately (READ 0) using a ViewLux uHTS microplate imager (PerkinElmer Life Sciences) equipped with excitation 525/20 and emission 598/25 filters. 2.5 μ l of a 2 \times Amplex Red solution (100 μ M Amplex Red (Cayman Chemical, Ann Arbor, MI), 200 μ M DTT (ThermoFisher Scientific), and 2 units/ml horseradish peroxidase (Sigma); diluted in HBSS and protected from light) were added to each well. Fluorescence was measured after a 15-min incubation at room temperature