

A systems biology approach combining ProLiFiler and Cancer Data Miner for an enhanced preclinical characterization of the WEE1 inhibitor adavosertib

Vincent Vuaroqueaux¹, Daniel Feger², Anne-Lise Peille¹, Oliver Siedentopf², Sadhana Panzade¹, Sarah Ulrich², Sebastian Dempe², Heinz-Herbert Fiebig¹, Jan Erik Ehlert²

¹4HF Biotec GmbH, Freiburg, Germany; ²Reaction Biology, Freiburg, Germany

Introduction

Adavosertib (AZD1775, MK-1775), is a clinical stage inhibitor of the tyrosine kinase Wee1. It plays an important role in regulating G2/M and S checkpoints. In case of DNA damage, Weel activation results in cell cycle arrest and mitosis entry failure until the DNA is repaired. In cancer cells, Wee1 inhibition leads to cell death per mitotic catastrophe.



In the present study, adavosertib was investigated in a cell survival and proliferation assay with 140 human cancer cell lines (CLs) representing all major tumor types (ProLiFiler platform of Reaction Biology) followed by mechanism of action (MoA) and biomarker analyses using 4HF Biotec's Cancer Data Mining in silico platform (DataMinerTM).

In vitro antitumor activity using ProLiFiler

> Assay procedure of the Cell Proliferation Assay

The ProLiFiler assay is performed using a contact-free nano-dispensing system (Tecan D300E) requiring small amounts of a test compound.



Assay procedure: Cells are seeded into 384 multi-well plates (A). The next day, compounds are added (B), and (C). Subsequently, the cell viability dve CellTiter-Glo luminescence is measured as a parameter for cell viability (D). IC₅₀ values are determined from 8 concentrations in duplicates.

Origin of cell lines

The panel currently consists of 140 human tumor cell lines derived from 21 tumor types.

Myelogenous Leukemia, DLBC: Lymphoma_Diffuse Large B Cell, ALCL: Anaplastic large cell lymp

Tumor type	#	Entities
Breast	13	4 ER ⁺ ; 5 ERBB2 ⁺ ; 4 TNBC
Ovary	9	
Uterus	6	2 cervix; 4 endometrium
Prostate	3	1 AR ⁺ ; 2 AR ⁻
Kidney	4	
Colorectal	11	
Stomach	9	
Pancreas	4	
Liver	2	2 hepatocellular carcinoma
Non-S mall Cell Lung	27	17 adenocarcinoma; 5 squamous; 3 large cell; 2 unclassified
Small Cell Lung	4	
Melanoma	4	
Nervous System	10	5 glioblastoma; 5 neuroblastoma
Sarcoma	6	3 osteosarcoma; 2 Ewing; 1 soft-tissue
Miscellaneous	4	1 bladder; 1 duodenum, 1 head & neck; 1 vulva (skin)
Leukemia	16	*3 ALL; 10 AML; 2 CLL; 1 CML
Myeloma	5	
Lymphoma	3	*2 DLBC; 1 ALCL
Total	140	
*ALL: Acute Lymphoblastic Leul	kemia, AML :	: Acute Myeloid Leukemia, CLL: Chronic Lymphocytic Leukemia ,CML: Chronic

(CLs: A-431, HUTU-80) Misce Sarcon
Le
Non Small Cell Lung ep
Uterus endo
Small Cell Lung
Liver hepat
M
Osteo
Sarcoma so
Co
F
Non Small Cell Lung La
Head a
Non Small Cell Lung adenocar
Non Small Cell Lung und
Neurol
Head & Neck, Stomach. (

🔲 Kidney & Bladder		L
Hematological		S
Brain		Ν
	_	

Drug Mechanism of Action (MoA)

The MoA Finder tool aims to identify the MoA of test compounds based on the results obtained in the ProLiFiler cell panel screening. To this end, we compare the sensitivity profile of a drug candidate to the profiles of more than 900 drugs for which the MoA is known.

- The MoA Finder tool allows the confirmation of the mechanism action of drugs.
- The data help to identify potential offtargets and to predict side effects.
- The MoA of compounds with unknown mechanisms, potential MoA can be predicted and suggestions for target deconvolution in a phenotypic approach can be made.

Database: The drug sensitivity profiles which reflect the MoAs of more than 900 anticancer agents, were collected from both the literature and/or determined in internal experiments.

Approach: The drug sensitivity profile (IC_{50}) from ProLiFiler screen) of a tes compound will be compared to the reterence compounds' drug sensitivit profiles. The correlation of the two datasets is evaluated by two statistical values, the Spearman Rho correlation value, and its pvalue.

Deliverables: Graphical presentation of the correlation analysis. Excel table with correlation results.

ProLiFiler – anti-proliferative activity of adavosertib



MoA Finder – drug mechanism of action of adavosertib

Non and

DNA repair

Apoptosi

Best correlating drug datasets

Inactive

cell cyle arrest



Biological context of Wee1 inhibition

The 10 Top-ranked MoA and correlation results (> 900 tested)*;**.

Rank	Compound	Mechanism of Action	Targets	₽ (Spearman test)	BH* adjusted p-value	Cell line (n=)	Geo. Mean (µM)	Database
1	Adavosertib	Cell cycle, genome integrity	<u>WEE1,</u> PLK1	0.57	7.76E-07	94	1.94	GDSC2
2	YK-4-279	DNA/RNA synthesis	EWS-FLI1, DHX9, ERG, ETV1	0.46	7.81E-04	95	4.65	GDSC1
3	AZD7762	Cell cycle, genome integrity	<u>СНЕК1</u> , СНЕК2	0.44	1.17E-03	94	1.07	GDSC2
4	Wee1 Inhibitor	Cell cycle genome integrity	<u>WEE1,</u> CHEK1	0.42	2.36E-03	93	8.63	GDSC2
5	AP-24534	ABL signaling	BCR-ABL, FGFR, PDGFR, VEGFR, SRC	0.4	2.13E-03	108	1.5	GDSC1
6	Pevonedistat	Cell cycle, genome integrity	NAE	0.4	5.76E-03	92	2.36	GDSC2
7	AZD6738	Cell cycle, genome integrity	ATR	0.39	6.79E-03	92	8.34	GDSC2
8	YK-4-279	DNA/RNA synthesis	EWS-FLI1, DHX9, ERG, ETV1	0.38	9.47E-03	92	10.92	GDSC2
9	Pevonedistat	Cell cycle, genome integrity	NAE	0.38	1.05E-02	86	1.24	GDSC1
10	VE821	Cell cycle, genome integrity	ATR/ATM	0.37	1.02E-02	90	57.94	GDSC2

One to one comparison with other MoAs



Figure 2. Top left: Volcano plot representation of MoA finder results of 471 anti-cancer drugs profiles (IC₅₀ data) with that of adavosertib, which is based on the IC₅₀ data from the ProLiFiler (COMPARE analysis*;**). Top right: Table showing the correlation results of the 10 top-ranked compounds. Bottom left: Graphical representation of the biological context of the proteins inhibited by TOP correlating drugs. Bottom right: Scatter plot showing the correlation between the sensitivity profiles of adavosertib and talazoparib (PARP inhibitor).*Huang, R., Wallqvist, A. & Covell, D.G. Anticancer metal compounds in NCI's tumor-screening database: putative mode of action. Biochem Pharmacol, (2005). Paull, K.D. et al. Display and analysis of patterns of differential activity of drugs against human tumor cell lines: development of mean graph and COMPARE algorithm. J Natl Cancer Inst, (1989). **Only drug datasets with more than 80 cell lines for comparison are shown in the table. 4HF Biotec (n=45), Sanger GDSC1 (n=251), Sanger GDSC2 (n=175).

* BH: Benjamini & Hochberg

Biomarker Analysis

The Biomarker Analysis tool reveals the specific genomic signature of tumor cell lines that are sensitive to a given test compound. To this end, we correlate the drug sensitivity profile (IC₅₀ data of ProLiFiler screen) of a test drug with available gene expression and genetic alterations (mutations, deletions, amplifications) datasets of the tumor cell lines.

The Biomarker Analysis tool can be used to identify genes expression changes or genetic alterations whose presence is indicative of the efficacy of the test compound and may serve as predictive biomarkers in disease models and patient cohorts

Approach: The ProLiFiler results are correlated with the gene expression, whole-exome sequencing (mutations), and somatic copy number alteration data of the tumor cell lines. An analysis is performed for every tumor cell line to determine the correlation of its drug sensitivity to each gene with regards to the expression level, mutation status, and somatic copy number.



Genomic alterations predicting sensitivity to adavosertib (using the GDSC2 dataset)

Mutations

Volcano plot showing gene mutations most significantly associated with tumor sensitivity to adavosertib (Wilcoxon test)



Amplifications

Volcano plot showing gene amplifications most significantly associated with tumor sensitivity to adavosertib (Wilcoxon test)



Testing set

	All cancers (Hematological & Solid)	Molecular signatures (MSigDB) containing genes highly expressed in	Molecular signatures (MSigDB) containing genes highly expressed in	Genes and pathways highly represented in CLs sensitive to adavosertib	Genes and pathways highly represented in CLs resistant to adavosertib
Dataset	ProLiFiler	CLs sensitive to adavosertib	CLs resistant to adavosertib	Train of all of the	A Station of
Cell lines with adavosertib data +gene expression profile (n=)	140	Myc Targets V1 *5.34e-13 E2F Targets *4.75e-08 Myc Targets V2 *5.79e-07 G2-M Checkpoint *1.71e-05	Estrogen Response Early *2.66e-14 Estrogen Response Late *6.80e-10 Epithelial Mesenchymal Transition *6.80e-10 Glycolysis *9.21e-09	NCM4 MCM5 MCM5 MCM5 MCM5 MCM5 MCM5 MCM5 M	RETSAT KRT18 CLIC3
Number of genes with Spearman rho >0 and p-values <0.05 (n=)	1141	DNA Repair *2.91e-03 Allograft Rejection *4.61e-03 IL-6/JAK/STAT3 Signaling 1.01e-02 Unfolded Protein Response 5.12e-02	Protein Secretion *7.832-08 UV Response Dn *5.49e-07 Apoptosis *1.38e-06 Xenobiotic Metabolism *2.68e-05	B POSB MINBAI MCMA2 NUDT21 POLD1 DDX18 EXXSC5 HCLS1	PROSESS CCND1 BLVRB PAPSS2 IGFBP4 ACOX2 FN1 SERPINE
Number of genes with S pearman rho <0 and p-values <0.05 (n=)	989	Wnt-beta Catenin Signaling 5.36e-02IL-2/STAT5 Signaling 6.61e-02024 $\begin{array}{c} 6\\ -\log 10 \\ (p-value) \end{array}$ 1012	Complement *4.49e-04 Coagulation *8.46e-04 0 2 4 6 8 10 12 14	IFNOR2 TGFB1 IRF8 GPR65 CCND2 PIM1	CFB CLU PLAT GPC1 NT6E PLOD2 RGS4

Validation sets

	Hematologi	S	
Dataset	GDSC2	CT R D ²	GD
Cell lines with adavosertib data +gene expression profile (n=)	88	137	4
Number of genes with Spearman rho >0 and p-values <0.05 (n=)	827	606	11
Number of genes with Spearman rho <0 and p-values <0.05 (n=)	1143	1413	22
	Shared significant genes (GDSC2 +CTD ²)		Shar ((
Number of genes with Spearman rho >0 and p-values <0.05 (n=)	184		
Number of genes with Spearman rho <0 and	3	14	

Figure 3: Left: Volcano plots showing levels of significance of individual genes for the association between adavosertib IC50 values and their gene mutation status (top plot) and gene amplification/deletions (bottom plot) (testing dataset: GDSC2; n = 761). The red and green dots show the gene alterations significantly associated with the response to adavosertib (Wilcoxon test; p-value < 0.05). Right: Transcripts correlating with CL response to adavosertib using ProliFiler dataset (top) along with validation in external datasets (bottom). Tables presenting the cell line panel used for the analysis for each dataset and the number of genes / probe sets significantly positively and negatively correlated with the adavosertib IC50 values (Spearman correlation test). The bar plots represent the results of the gene set enrichment analysis (Molecular Signatures Database Hallmark 2020) of genes correlating with adavosertib IC50 values in ProfiLer and external datasets. Clustergrams show the genes underlying in the 10 top-ranked gene expression signatures of the MSigDB Hallmarks

Xie Z, et al., Gene set knowledge discovery with Enrichr. Current Protocols (2021). Liberzon A et al., The Molecular Signatures Database Hallmark Gene Set Collection. Cell Systems (2015).

In conclusion, our studies demonstrate broad anticancer activity of adavosertib and confirm its proposed mechanism of action. The biomarkers we identified will facilitate the selection of pre-clinical in vivo tumor models and, if confirmed, even patient selection for clinical trials. The combined use of the ProLiFiler and Cancer Data Miner Platforms has the potential to accelerate and de-risk the development of anti-cancer agents.



Biomarker Screening

Transcripts and gene signatures predicting cancer sensitivity to adavosertib



Conclusions