

BRD2 Bromodomain Assay Service (ThermoShift)

BRD2-Tndm-His [Reader Domain Assay](#) for compound screening and profiling using the thermoshift assay format to measure compound binding to the target.

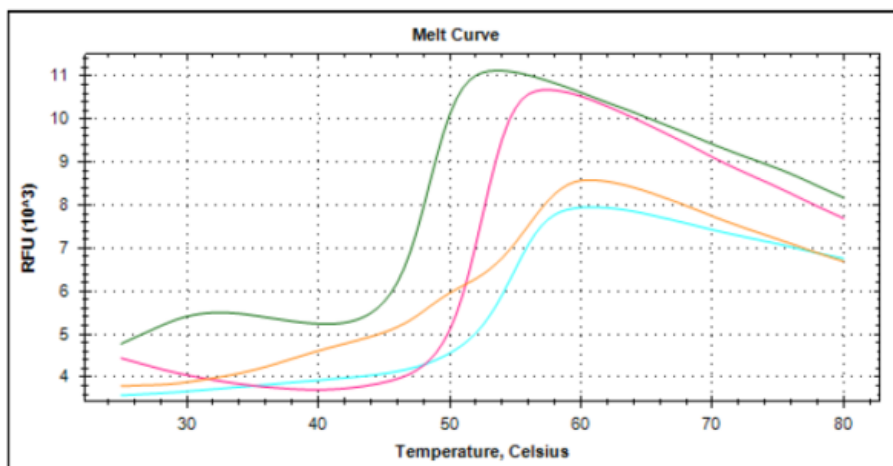
Target Overview

CAT#	Tm-BRD2-T-H
Protein	Human BRD2-T bromodomain, His-tagged
Synonyms	O27.1.1, RING3

Assay Properties

Readout	Thermal stability change after ligand binding
Concentration of protein in assay	0.125 mg/ml (2.3 μ M)
Concentration of compound in assay	25 μ M
Apo form T_m ($^{\circ}$C)	48,5
Reference Compound T_m s:	JQ1 - 7.5 (Orange)PFI1 - 4 (Pink)BET151 - 6.5 (Blue)No stabilization was observed in the presence of other control compounds (CBP112, RVX-208, GSK2801, and PFI3)
Scientific Information	P25440 (http://www.uniprot.org/uniprot/P25440)
Screening Location	Malvern, USA
Further information	More information can be found on our website Reader Domain Assay Services .

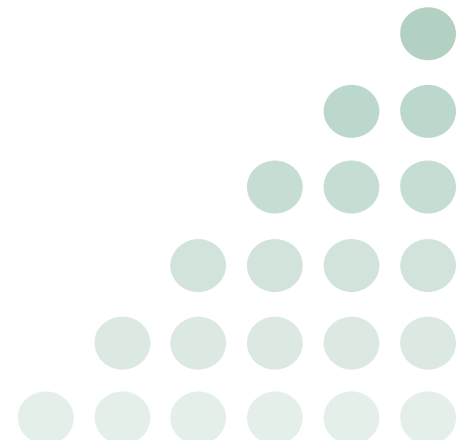
Melt Curve



Let's Discover Together.

© Oct 2022 **Reaction Biology**. All rights reserved.

https://www.reactionbiology.com/datasheet/brd2-tndm-hisreader_thermoshift_malvern



Melt Peak

