

BRDT-1 Bromodomain Assay Service (ThermoShift)

BRDT-1-His [Reader Domain Assay](#) for compound screening and profiling using the thermoshift assay format to measure compound binding to the target.

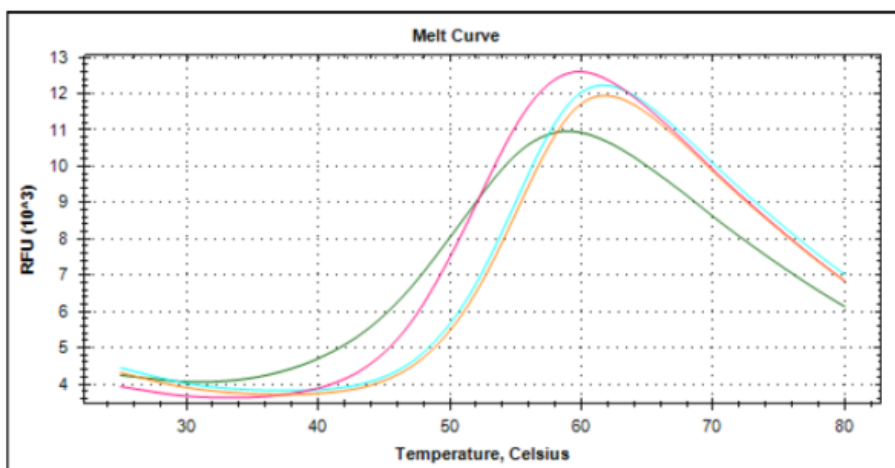
Target Overview

CAT#	Tm-BRDT-1-H
Protein	Human BRDT-1 bromodomain, His-tagged
Synonyms	CT9, RING3-like protein, BRD6

Assay Properties

Readout	Thermal stability change after ligand binding
Concentration of protein in assay	0.075 mg/ml (4.5 μ M)
Concentration of compound in assay	25 μ M
Apo form T_m ($^{\circ}$C)	51,5
Reference Compound T_m s:	JQ1 - 3.5 (Orange)PFI1 - 1 (Pink)BET151 - 3.5 (Blue)No stabilization was observed in the presence of other control compounds (CBP112, SGC-CBP30, RVX-208, GSK2801, and PFI3)
Scientific Information	Q58F21 (http://www.uniprot.org/uniprot/Q58F21)
Screening Location	Malvern, USA
Further information	More information can be found on our website Reader Domain Assay Services .

Melt Curve



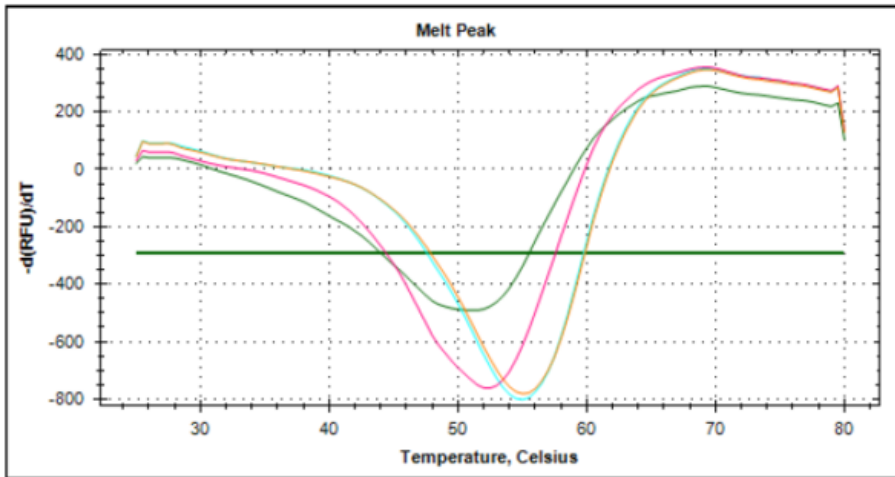
Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/brdt-1-hisreader_thermoshift_malvern



Melt Peak



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/brdt-1-hisreader_thermoshift_malvern

