

BRD3-2 Bromodomain Assay Service (ThermoShift)

BRD3-2-His [Reader Domain Assay](#) for compound screening and profiling using the thermoshift assay format to measure compound binding to the target.

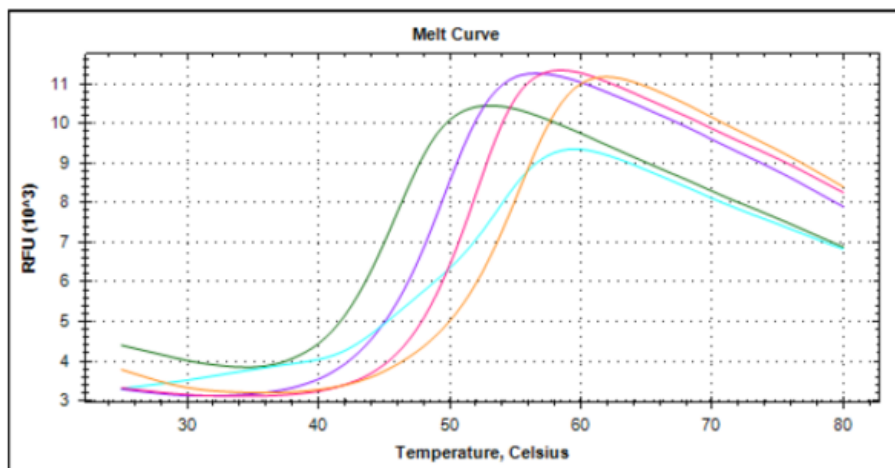
Target Overview

CAT#	Tm-BRD3-2-H
Protein	Human BRD3-2 bromodomain, His-tagged
Synonyms	RING3-like protein

Assay Properties

Readout	Thermal stability change after ligand binding
Concentration of protein in assay	0.085 mg/ml (5.4 μ M)
Concentration of compound in assay	25 μ M
Apo form T_m ($^{\circ}$ C)	46
Reference Compound T_m s:	JQ1 - 9.5 (Orange)PFI1 - 6 (Pink)BET151 - 7.5 (Blue)RVX-208 - 3.5 (Purple)
Scientific Information	Q15059 (http://www.uniprot.org/uniprot/Q15059)
Screening Location	Malvern, USA
Further information	More information can be found on our website Reader Domain Assay Services .

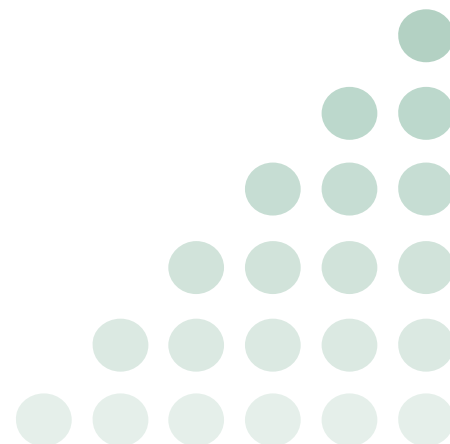
Melt Curve



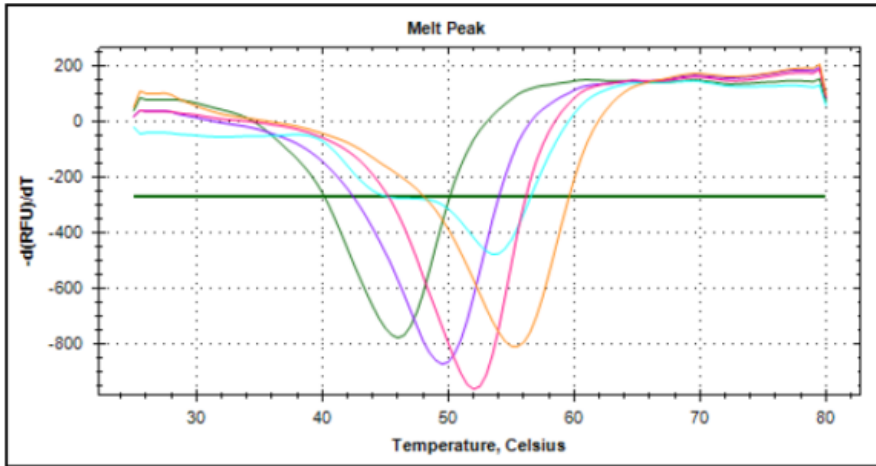
Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/brd3-2-hisreader_thermoshift_malvern



Melt Peak



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/brd3-2-hisreader_thermoshift_malvern

