

BRD1 Bromodomain Assay Service (ThermoShift)

BRD1-GST [Reader Domain Assay](#) for compound screening and profiling using the thermoshift assay format to measure compound binding to the target.

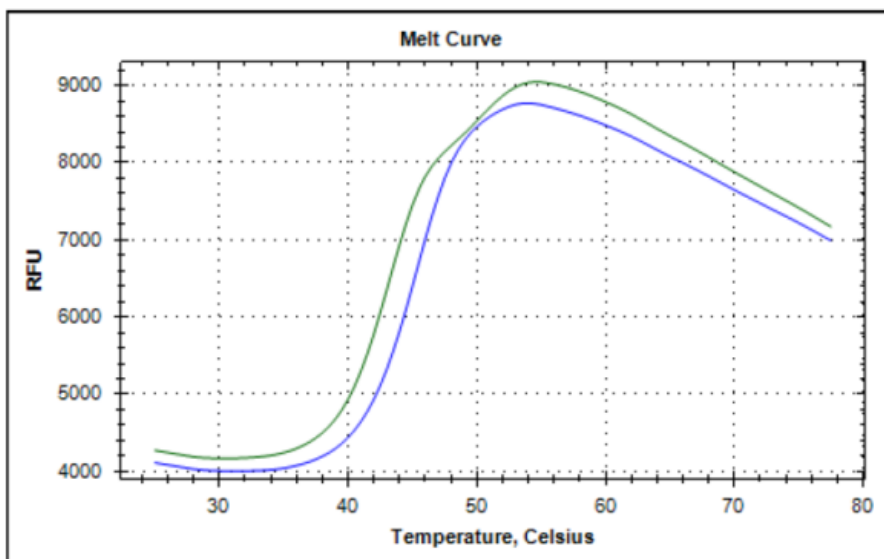
Target Overview

CAT#	Tm-BRD1-GST
Protein	Human BRD1 bromodomain, GST-tagged
Synonyms	BR140-like protein, BRL; Bromodomain and PHD finger-containing protein 2, BRPF-2

Assay Properties

Readout	Thermal stability change after ligand binding
Concentration of protein in assay	0.15 mg/ml (3.53 μ M)
Concentration of compound in assay	25 μ M
Apo form T_m ($^{\circ}$C)	43
Reference Compound T_m s:	Bromosporine - 2.5 (Blue) No stabilization was observed in the presence of other control compounds (JQ1, PFI1, CBP112, SGC-CBP30, BET151, RVX-208, GSK2801 and PFI3)
Scientific Information	O95696 (http://www.uniprot.org/uniprot/O95696)
Screening Location	Malvern, USA
Further information	More information can be found on our website Reader Domain Assay Services .

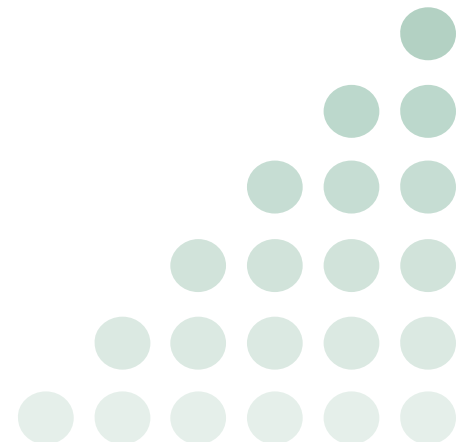
Melt Curve



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/brd1-greader_thermoshift_malvern



Melt Peak

