

BRWD1-2 Bromodomain Assay Service (ThermoShift)

BRWD1-2-GST [Reader Domain Assay](#) for compound screening and profiling using the thermoshift assay format to measure compound binding to the target.

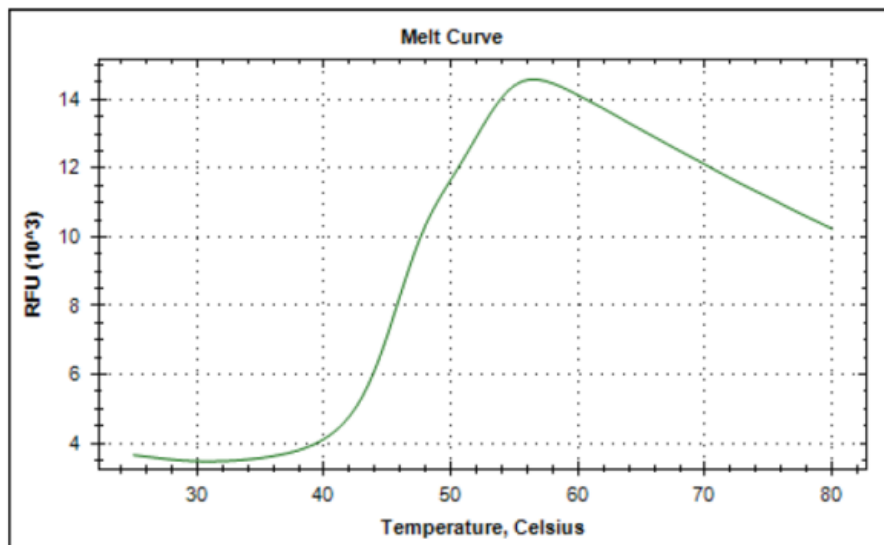
Target Overview

CAT#	Tm-BRWD1-2-GST
Protein	Human BRWD1-2 bromodomain, GST-tagged
Synonyms	Bromodomain and WD repeat-containing protein 1; WDR9

Assay Properties

Readout	Thermal stability change after ligand binding
Concentration of protein in assay	0.1 mg/ml (2.4 μ M)
Concentration of compound in assay	25 μ M
Apo form T_m ($^{\circ}$C)	46
Reference Compound T_m s:	No stabilization was observed in the presence of control compounds (JQ1, PFI1, CBP112, Bromosporine, SGC-CBP30, BET151, RVX-208, GSK2801 and PFI3)
Scientific Information	Q9NSI6 (http://www.uniprot.org/uniprot/Q9NSI6)
Screening Location	Malvern, USA
Further information	More information can be found on our website Reader Domain Assay Services .

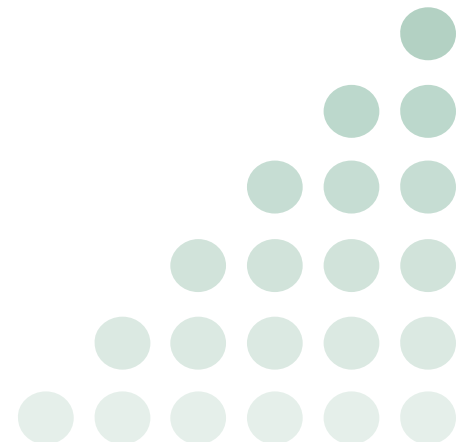
Melt Curve



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/brwd1-2-gstreader_thermoshift_malvern



Melt Peak

