

SP100 Bromodomain Assay Service (ThermoShift)

SP100-GST [Reader Domain Assay](#) for compound screening and profiling using the thermoshift assay format to measure compound binding to the target.

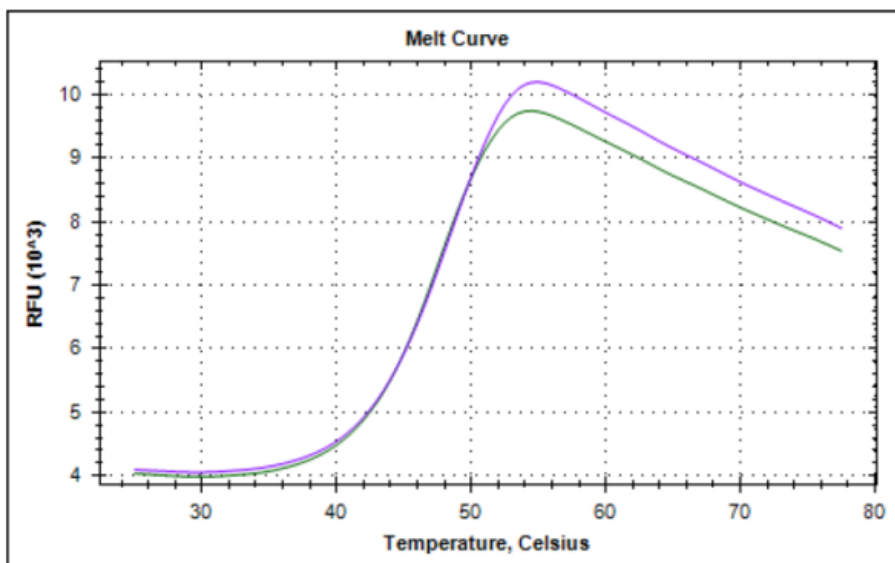
Target Overview

CAT#	Tm-SP100-GST
Protein	Human SP100 bromodomain, GST-tagged
Synonyms	Nuclear dot-associated Sp100 protein

Assay Properties

Readout	Thermal stability change after ligand binding
Concentration of protein in assay	0.1 mg/ml (2.08 μ M)
Concentration of compound in assay	25 μ M
Apo form T_m ($^{\circ}$C)	47,5
Reference Compound T_m s:	RVX-208 - 1 (Purple) No stabilization was observed in the presence of other control compounds (JQ1, PFI1, CBP112, Bromosporine, SGC-CBP30, BET151, GSK2801 and PFI3)
Scientific Information	P23497 (http://www.uniprot.org/uniprot/P23497)
Screening Location	Malvern, USA
Further information	More information can be found on our website Reader Domain Assay Services .

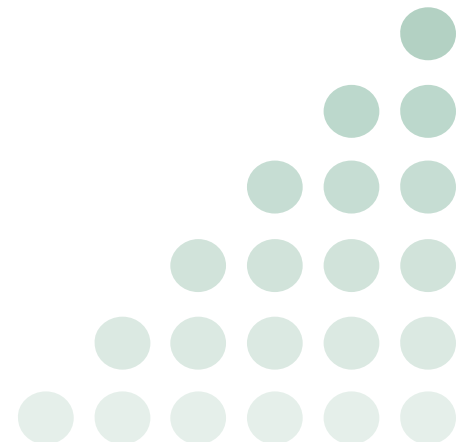
Melt Curve



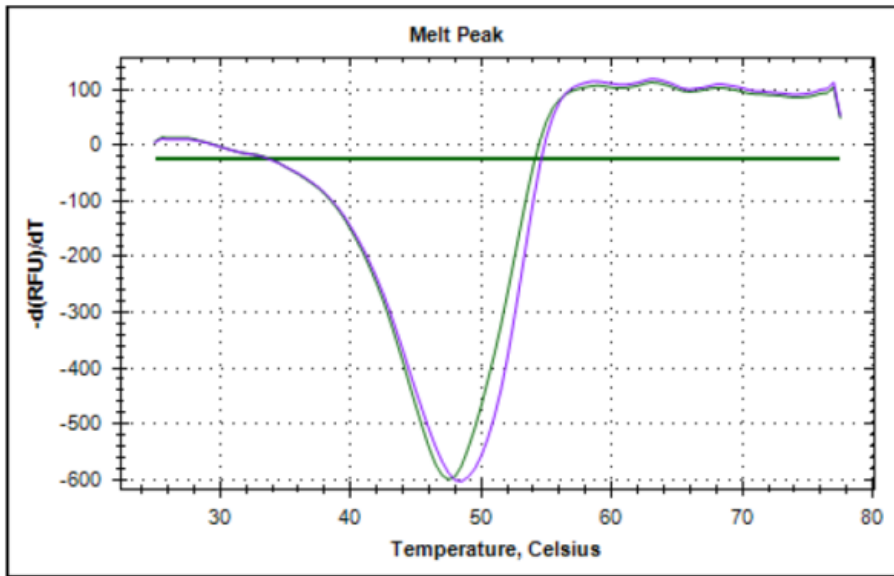
Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/sp100-gstreader_thermoshift_malvern



Melt Peak



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/sp100-gstreader_thermoshift_malvern

