

BAZ2B Bromodomain Assay Service (ThermoShift)

BAZ2B-His [Reader Domain Assay](#) for compound screening and profiling using the thermoshift assay format to measure compound binding to the target.

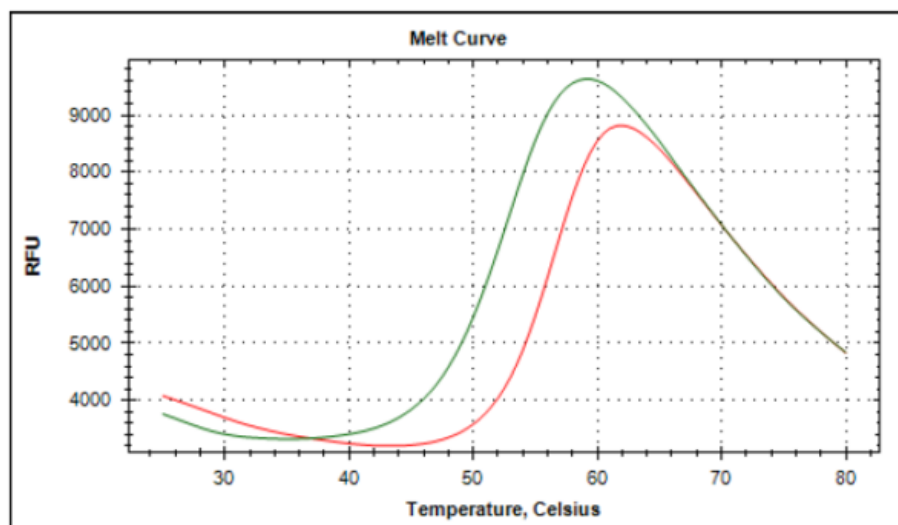
Target Overview

CAT#	Tm-BAZ2B-H
Protein	Human BAZ2B bromodomain, his-tagged
Synonyms	Bromodomain adjacent to zinc finger domain protein 2B; hWALp4; KIAA1476

Assay Properties

Readout	Thermal stability change after ligand binding
Concentration of protein in assay	0.1 mg/ml (6.2 μ M)
Concentration of compound in assay	25 μ M
Apo form T_m ($^{\circ}$C)	53
Reference Compound T_m s:	GSK2801 - 3.5 (red) No stabilization was observed in the presence of additional control compounds (JQ1, PFI1, CBP112, Bromosporine, SGC-CBP30, BET151, RVX-208 and PFI3)
Scientific Information	Q9UIF8 (http://www.uniprot.org/uniprot/Q9UIF8)
Screening Location	Malvern, USA
Further information	More information can be found on our website Reader Domain Assay Services .

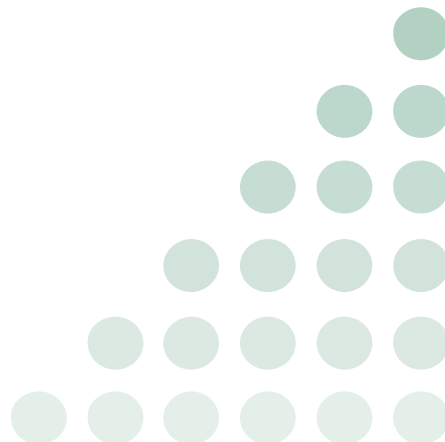
Melt Curve



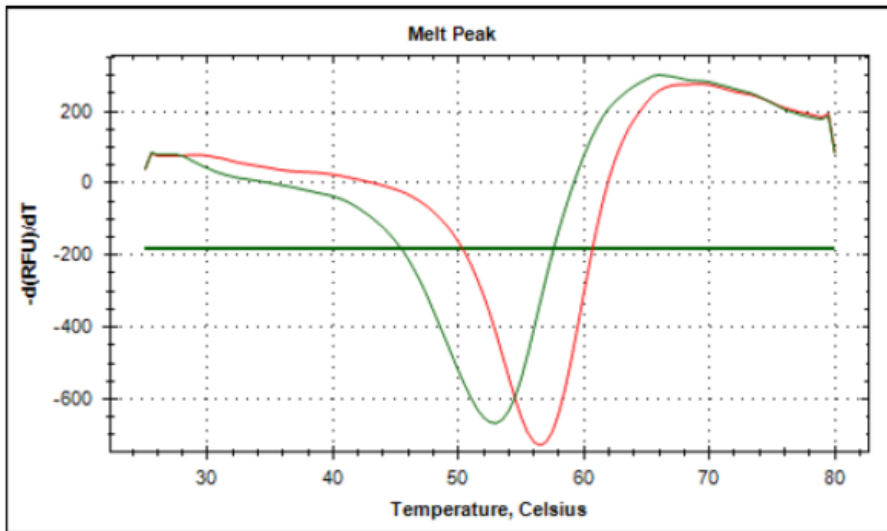
Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/baz2b-hreader_thermoshift_malvern



Melt Peak



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/baz2b-hreader_thermoshift_malvern

