

TRIM33b Bromodomain Assay Service (ThermoShift)

TRIM33b-His [Reader Domain Assay](#) for compound screening and profiling using the thermoshift assay format to measure compound binding to the target.

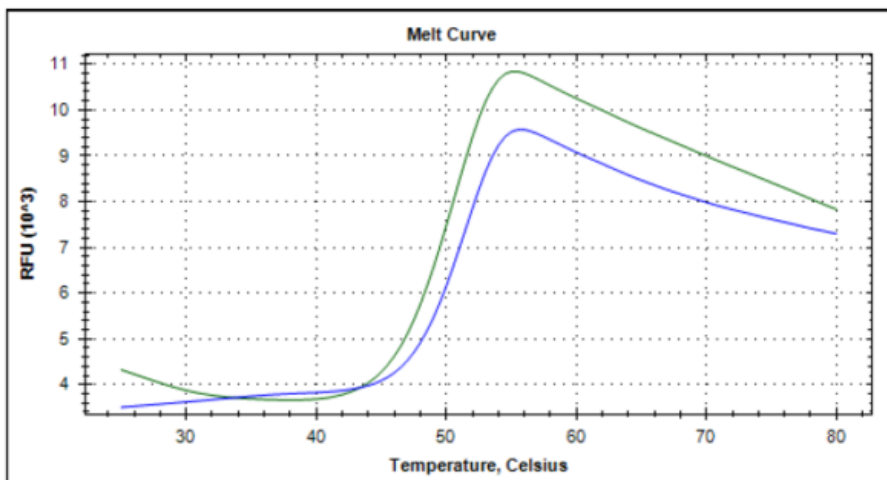
Target Overview

CAT#	Tm-TRIM33b-H
Protein	Human TRIM33b bromodomain, His-tagged
Synonyms	E3 ubiquitin-protein ligase TRIM33; Ectodermin homolog; Protein Rfg7; TIF1-gamma; Tripartite motif-containing protein 33

Assay Properties

Readout	Thermal stability change after ligand binding
Concentration of protein in assay	0.075 mg/ml (3.1 μ M)
Concentration of compound in assay	25 μ M
Apo form T_m ($^{\circ}$C)	50,5
Reference Compound T_m s:	Bromosporine - 1.0 (Blue) No stabilization was observed in the presence of other control compounds (JQ1, PFI1, CBP112, SGC-CBP30, BET151, GSK2801 and PFI3)
Scientific Information	Q9UPN9 (http://www.uniprot.org/uniprot/Q9UPN9)
Screening Location	Malvern, USA
Further information	More information can be found on our website Reader Domain Assay Services .

Melt Curve



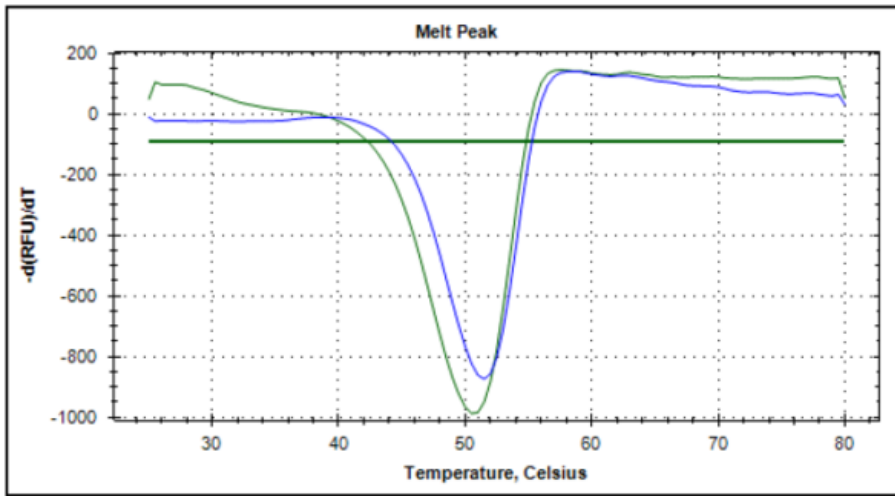
Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/trim33b-hreader_thermoshift_malvern



Melt Peak



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/trim33b-hreader_thermoshift_malvern

