

# BRPF3 Bromodomain Assay Service (ThermoShift)

BRPF3-GST [Reader Domain Assay](#) for compound screening and profiling using the thermoshift assay format to measure compound binding to the target.

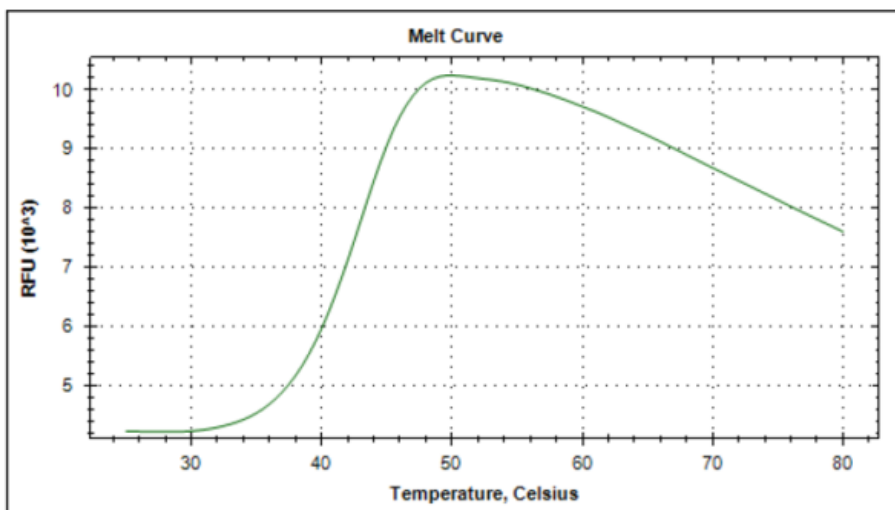
## Target Overview

CAT#	Tm-BRPF3-GST
Protein	Human BRPF3 bromodomain, GST-tagged
Synonyms	KIAA1286

## Assay Properties

Readout	Thermal stability change after ligand binding
Concentration of protein in assay	0.15 mg/ml (3.69 $\mu$ M)
Concentration of compound in assay	25 $\mu$ M
Apo form $T_m$ ( $^{\circ}$ C)	43
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	<a href="http://www.uniprot.org/uniprot/Q9ULD4">Q9ULD4</a> ( <a href="http://www.uniprot.org/uniprot/Q9ULD4">http://www.uniprot.org/uniprot/Q9ULD4</a> )
Screening Location	Malvern, USA
Further information	More information can be found on our website <a href="#">Reader Domain Assay Services</a> .

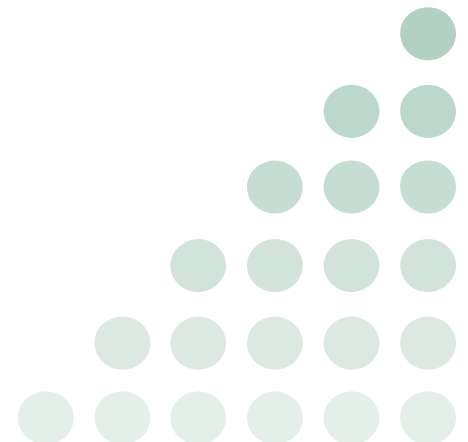
## Melt Curve



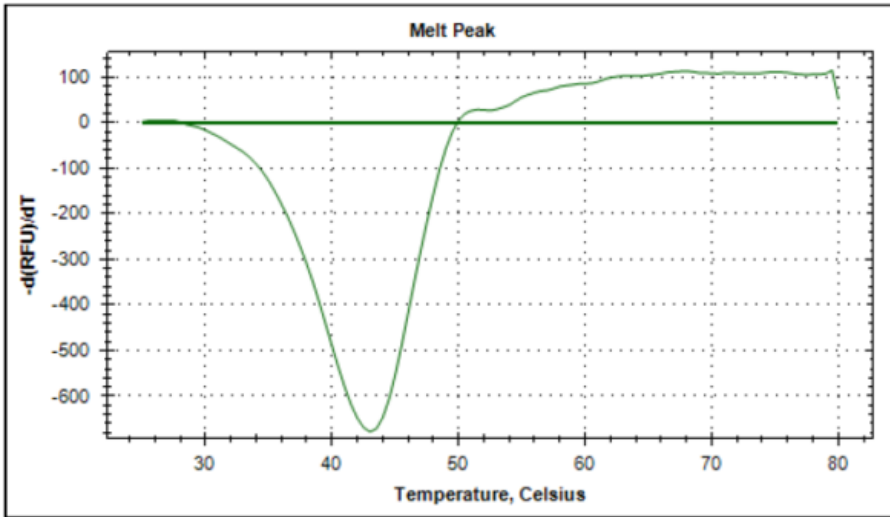
Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/brpf3-gstreader\\_thermoshift\\_malvern](https://www.reactionbiology.com/datasheet/brpf3-gstreader_thermoshift_malvern)



Melt Peak



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/brpf3-gstreader\\_thermoshift\\_malvern](https://www.reactionbiology.com/datasheet/brpf3-gstreader_thermoshift_malvern)

