

CREBBP Bromodomain Assay Service (ThermoShift)

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

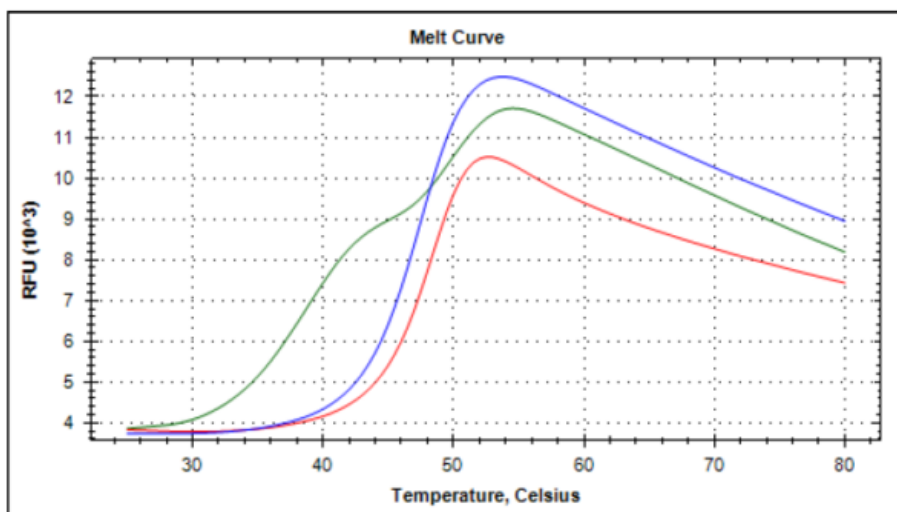
Target Overview

CAT#	Tm-CREBBP-GST
Protein	Human CREBBP bromodomain, GST-tagged
Synonyms	CREB-binding protein, CBP

Assay Properties

Readout	Thermal stability change after ligand binding
Concentration of protein in assay	0.1 mg/ml (2.44 μ M)
Concentration of compound in assay	25 μ M
Apo form T_m ($^{\circ}$C)	39
Reference Compound T_m s:	CBP112 - 8.5 (Blue)SGC-CBP30 - 9.5 (Red)No stabilization was observed in the presence of other control compounds (JQ1, RVX-208, GSK2801 and PFI3)
Scientific Information	Q92793 (http://www.uniprot.org/uniprot/Q92793)
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/crebbp-gstreader_thermoshift_malvern



Melt Peak

