

SMARCA4 Bromodomain Assay Service (ThermoShift)

SMARCA4-His [Reader Domain Assay](#) for compound screening and profiling using the thermoshift assay format to measure compound binding to the target.

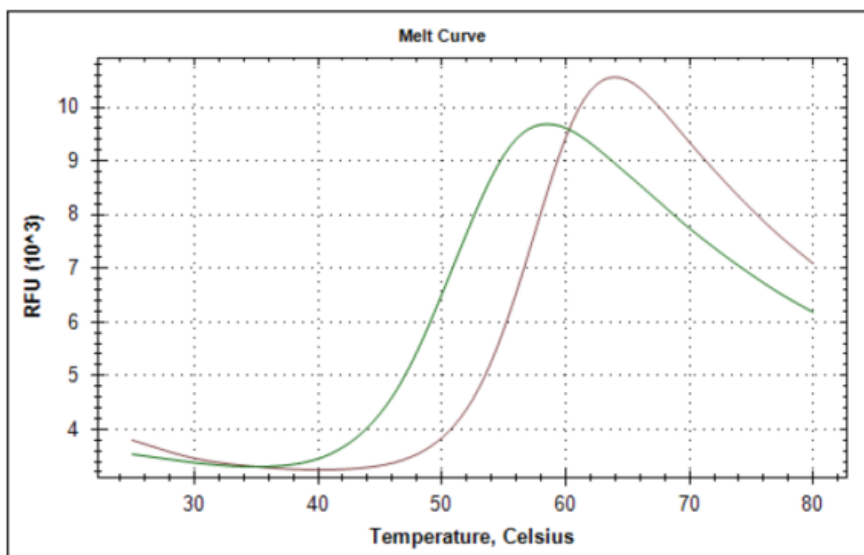
Target Overview

CAT#	Tm-SMARCA4-H
Protein	Human SMARCA4 bromodomain, His-tagged
Synonyms	BRG1, BAF190A, Protein BRG-1, SNF2-beta

Assay Properties

Readout	Thermal stability change after ligand binding
Concentration of protein in assay	0.15 mg/ml (8.8 μ M)
Concentration of compound in assay	25 μ M
Apo form T_m ($^{\circ}$C)	51,5
Reference Compound T_m s:	PFI3 - 6 (Purple) No stabilization was observed in the presence of other control compounds (JQ1, PFI1, CBP112, Bromosporine, SGC-CBP30, BET151, RVX-208, and GSK2801)
Scientific Information	P51532 (http://www.uniprot.org/uniprot/P51532)
Screening Location	Malvern, USA
Further information	More information can be found on our website Reader Domain Assay Services .

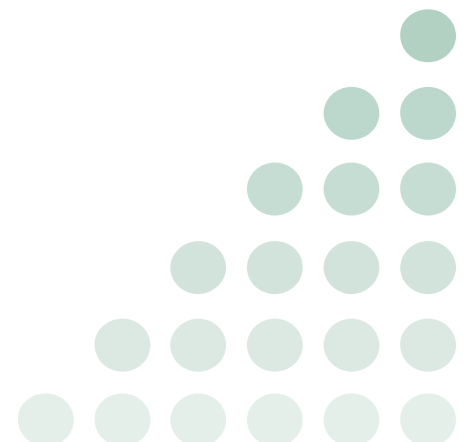
Melt Curve



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/smarca4-hisreader_thermoshift_malvern



Melt Peak

