

SP140 Bromodomain Assay Service (ThermoShift)

SP140-GST [Reader Domain Assay](#) for compound screening and profiling using the thermoshift assay format to measure compound binding to the target.

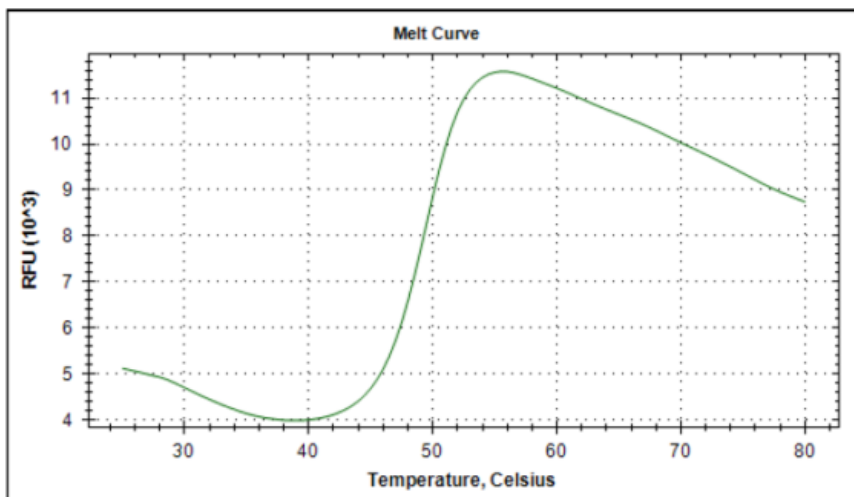
Target Overview

| | |
|-----------------|--|
| CAT# | Tm-SP140-GST |
| Protein | Human SP140 bromodomain, GST-tagged |
| Synonyms | Lymphoid-restricted homolog of Sp100; Nuclear autoantigen Sp-140; Speckled 140 kDa |

Assay Properties

| | |
|---|--|
| Readout | Thermal stability change after ligand binding |
| Concentration of protein in assay | 0.1 mg/ml (2.1 μ M) |
| Concentration of compound in assay | 25 μ M |
| Apo form T_m ($^{\circ}$C) | 49,5 |
| 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 | Q13342 (http://www.uniprot.org/uniprot/Q13342) |
| 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/sp140-gstreader_thermoshift_malvern



Melt Peak

