

UHRF1-[TDR] (His) Reader Domain Assay Service

UHRF1-[TDR] (His) [Reader Domain Assay](#) for compound screening and profiling using the thermoshift assay format to measure compound binding to the target.

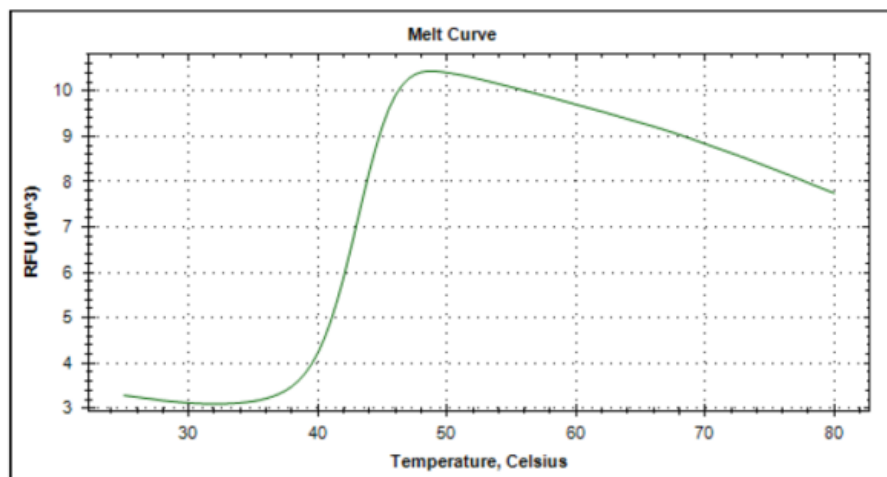
Target Overview

CAT#	Tm-UHRF1-[TDR]-H
Protein	Human UHRF1 tandem Tudor domain, his tagged
Synonyms	E3 ubiquitin-protein ligase UHRF1, Nuclear protein 95

Assay Properties

Readout	Thermal stability change after ligand binding
Concentration of protein in assay	0.1 mg/ml (4.5 μ M)
Concentration of compound in assay	N/A
Apo form T_m ($^{\circ}$C)	43
Reference Compound T_m s:	No reference compound is currently available for this protein
Scientific Information	Q96T88 (http://www.uniprot.org/uniprot/Q96T88)
Screening Location	Malvern, USA
Further information	More information can be found on our website Reader Domain Assay Services .

Melt Curve



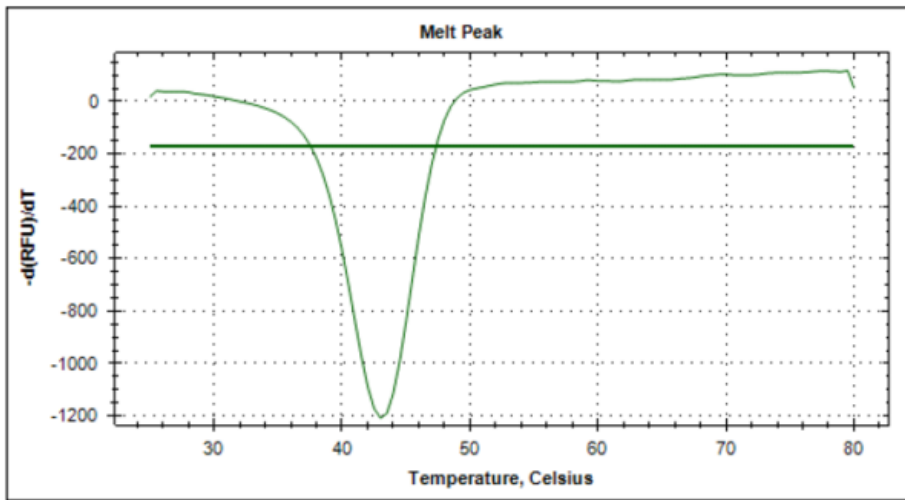
Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/uhrf1-tdr-hreader_thermoshift_malvern



Melt Peak



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/uhrf1-tdr-hreader_thermoshift_malvern

