

## UHRF1 [TDR-PHD] (His)

**CATALOG NO.:** RD-11-248

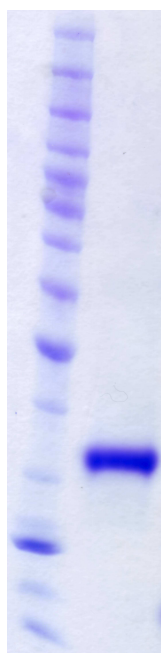
**LOT NO.:**

**DESCRIPTION:** Human recombinant UHRF1 [TDR-PHD] bromodomain (residues 123-366; Genbank Accession # NM\_001048201; MW = 31.0 kDa) expressed as an N-terminal His-fusion protein in *E. coli*.

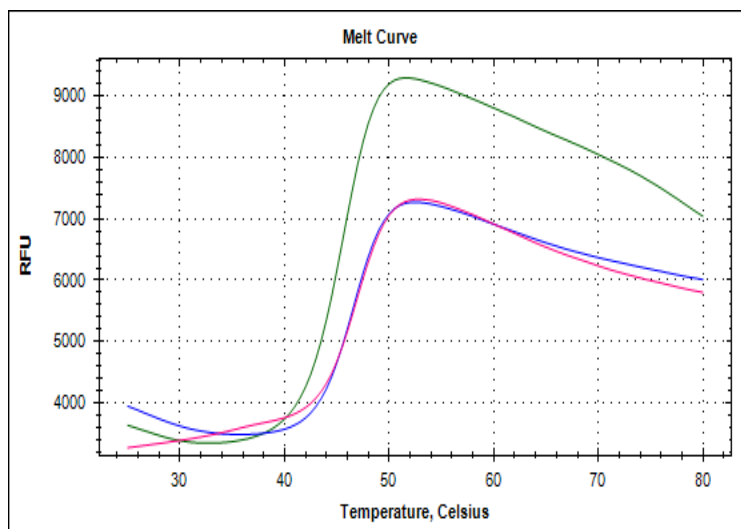
**PURITY:** >95% by SDS-PAGE

**SUPPLIED AS:**  $\_ \mu\text{g}/\mu\text{L}$  in 50 mM Tris HCl, pH 7.5, 500 mM NaCl, 1 mM TCEP, 10 % glycerol

**STORAGE:** -70°C. Thaw quickly and store on ice before use. The remaining, unused, undiluted protein should be snap frozen, for example in a dry/ice ethanol bath or liquid nitrogen. Minimize freeze/thaws if possible, but very low volume aliquots (<5  $\mu\text{l}$ ) or storage of diluted enzyme is not recommended.



**Coomassie blue-stained SDS-PAGE (12% acrylamide) of 5  $\mu\text{g}$  of RBC UHRF1-[TDR-PHD](His).** MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, **50**, 40, 30, 25, **20**, 15, 10 kDa.



**Differential Scanning Fluorimetry of RBC UHRF1[TDR-PHD](His)**  
Thermal denaturation of UHRF1[TDR-PHD](His) is detected (CFX384™ Touch thermal cycler, 'FRET' channel; Bio-Rad) by increased binding and fluorescence of the dye SYPRO® Orange (Life Technologies). Apo form of UHRF1[TDR-PHD](His) displays a  $T_m$  of 45.5°C. Addition of 25  $\mu\text{M}$  Bromosporine (blue) or PF11 (pink) stabilizes the protein folding and shifts the  $T_m$  (inflection point to 46.5°C or 47°C respectively).

This product is not intended for therapeutic or diagnostic use in animals or in humans.

Reaction Biology

1 Great Valley Parkway, Malvern PA, USA 19355

requests@reactionbiology.com www.reactionbiology.com