

UHRF1-[TDR-PHD] (GST)

CATALOG NO.: RD-11-300

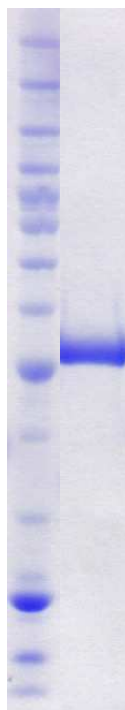
LOT NO.:

DESCRIPTION: Human recombinant UHRF1-[TDR-PHD](GST) bromodomain (residues 123-366; Genbank Accession # NM_001048201; MW = 55.3 kDa) expressed as an N-terminal GST-fusion protein in *E. coli*.

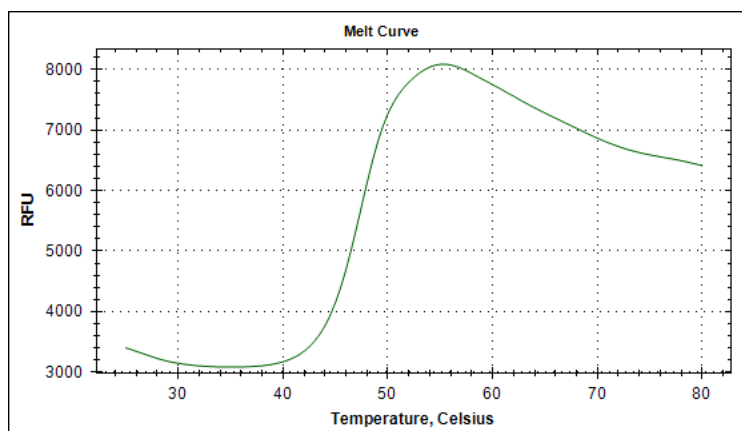
PURITY: >95% by SDS-PAGE

SUPPLIED AS: \sim μ g/ μ 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 μ L) or storage of diluted enzyme is not recommended.



Coomassie blue-stained SDS-PAGE (12% acrylamide) of 5 μ g of RBC UHRF1-[TDR-PHD] (GST). 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 Thermal denaturation of UHRF1-[TDR-PHD] (GST) is detected (CFX384TM 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]-(GST) displays a T_m of 47.5°C, and is not stabilized in the presence of various known bromodomain ligands (JQ1, PFI1, CBP112, Bromosporine, SGC-CBP30, BET151, RVX-208, GSK2801 and PFI3; all tested at 25 μ M).

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