

## UHRF1-[SRA] (His)

**CATALOG NO.:** RD-11-235

**LOT NO.:**

**DESCRIPTION:** Human recombinant UHRF1 SRA domain (residues 414-617; Genbank Accession # NM\_001048201; MW = 25.4 kDa) expressed as an N-terminal His-fusion protein in *E. coli*. A tumor promoter overexpressed in a wide variety of cancers, the epigenetic regulator UHRF1 (Ubiquitin-like PHD and Ring Finger 1) is a multidomain protein that co-localizes with DNMT1 and functions in the maintenance of DNA methylation patterns (see review<sup>1</sup>). Its SRA (SET and Ring-Associated) domain is an epigenetic reader that specifically binds 5-methylcytosine at hemimethylated CpG sites<sup>2-5</sup>.

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

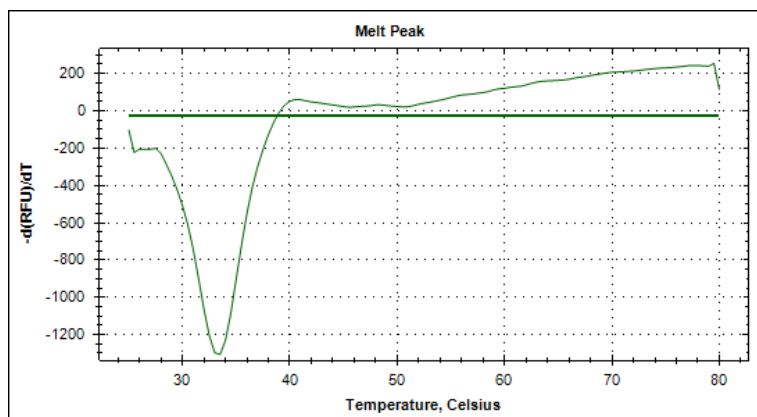
**SUPPLIED AS:**  $\_ \mu\text{g}/\mu\text{L}$  in 50 mM Tris HCl, pH 7.5, 500 mM NaCl, 1 mM TCEP, 10 % glycerol as determined by OD<sub>280</sub>.

**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.

**REFERENCES:** 1) W. Ashraf *et al. Oncotarget* 2017 **8** 51946; 2) L.M. Johnson *et al. Curr. Biol.* 2007 **17** 322; 3) G.V Avvkumov *et al. Nature* 2008 **455** 822; 4) C. Qian *et al. J. Biol. Chem.* 2008 **283** 34490; 5) J. Song & G.P. Pfeifer *Bioessays* 2016 **38** 1038



**Coomassie blue-stained SDS-PAGE (4-12% acrylamide) of 4  $\mu\text{g}$  of RBC UHRF1-[SRA] (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-[SRA] (His)** Thermal denaturation of UHRF1[SRA](His) is detected (CFX384™ Touch thermal cycler, 'FRET' channel; Bio-Rad) by increased binding and fluorescence of the dye SYPRO® Orange (Life Technologies). In this negative differential plot of its melt curve UHRF1-[SRA](His) the downward pointing 'peak' indicates a T<sub>m</sub> of 33.5°C.

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