

Recombinant UHRF1 protein

Catalog No: 81015

Lot No: 14417001

Expressed In: *E. coli*

Quantity: 50, 1000 µg

Concentration: 1 µg/µl

Source: Human

Buffer Contents: Recombinant UHRF1 protein is supplied at a concentration of 1 µg/µl in 25 mM Tris pH 8.0, 300 mM NaCl, 5% glycerol.

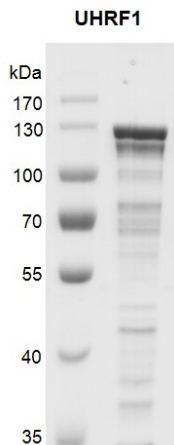
Background: UHRF1 (Ubiquitin Like With PHD And Ring Finger Domains 1, also known as Np95, hNP95, ICBP90, RNF106, TDRD22, hUHRF1, huNp95) is a member of a subfamily of RING-finger type E3 ubiquitin ligases. UHRF1 binds to specific DNA sequences, and recruits a histone deacetylase to regulate gene expression. UHRF1 plays a major role in the G1/S transition by regulating topoisomerase II α ; and retinoblastoma gene expression, and functions in the p53-dependent DNA damage checkpoint. UHRF1 is overexpressed in various cancers. UHRF1 plays a critical role in maintaining DNA methylation in mammalian cells.

Protein Details: Recombinant UHRF1 protein was expressed in *E. coli* cells as the full length protein (accession number NP_037414.3) with a C-terminal 6×His tag. The molecular weight of the protein is 93.1 kDa.

Application Notes:

This protein is useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.

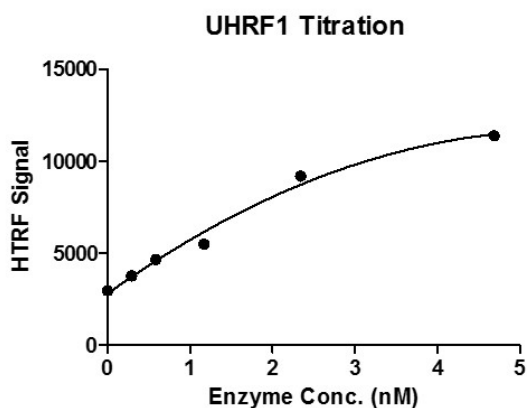


Recombinant UHRF1 protein gel

8% SDS-PAGE Coomassie staining

MW: 93.1 kDa

Purity: > 60%



UHRF1 binding assay 1 μ M H3K9me3 peptide was incubated with different concentrations of UHRF1 protein in 50 mM HEPES-NaOH pH 7.5, 0.1% BSA for 1 hour, then 10 μ l anti-His antibody and SA-XL665 mixture (1:100 dilution in the same buffer) was added to each reaction system and incubated for 30 min. All the operations and reactions were performed at room temperature. HTRF assay was used for detection.

