
UHRF1 Polyclonal Antibody

(Catalog #A-2714)

Background

UHRF1, also named as ICBP90, NP95 and RNF106, is a putative E3 ubiquitin-protein ligase. It may participate in methylation-dependent transcriptional regulation. UHRF1 binds to inverted 5'-CCAAT-3' box 2 in the TOP2A promoter, and activates TOP2A expression. It is important for G1/S transition and may be involved in DNA repair and chromosomal stability. UHRF1's ability to repress its direct target gene expression is dependent on PHD(UHRF1) binding to unmodified H3R2.

Description

UHRF1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q96T88

Purification

Affinity Purified

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 1-260 of human UHRF1 (NP_001041666.1)

Storage

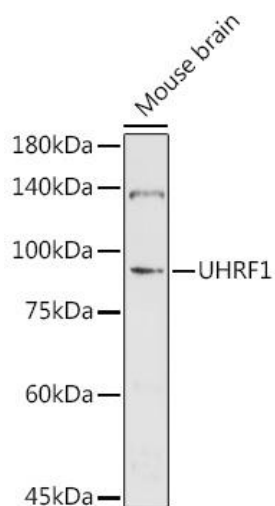
Shipped at 4°C. Upon receipt, store at -20°C. Avoid repeated freeze.

Alternative Names

FLJ21925, ICBP90, MGC138707, Np95, RNF106, hNP95

Application

ELISA, WB; Recommended dilution: WB 1:100 – 1:500



Western blot analysis of lysates from Mouse brain, using UHRF1 Polyclonal Antibody at 1:500 dilution.

Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.

Lysates/proteins: 25µg per lane.

Blocking buffer: 3% nonfat dry milk in TBST.

Exposure time: 90s.