
HUS1 Polyclonal Antibody

(Catalog # A71418)

Background

The protein encoded by this gene is a component of an evolutionarily conserved, genotoxin-activated checkpoint complex that is involved in the cell cycle arrest in response to DNA damage. This protein forms a heterotrimeric complex with checkpoint proteins RAD9 and RAD1. In response to DNA damage, the trimeric complex interacts with another protein complex consisting of checkpoint protein RAD17 and four small subunits of the replication factor C (RFC), which loads the combined complex onto the chromatin. The DNA damage induced chromatin binding has been shown to depend on the activation of the checkpoint kinase ATM, and is thought to be an early checkpoint signaling event. Alternative splicing results in multiple transcript variants.

Description

HUS1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

O60921

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 1-280 of human HUS1 (NP_004498.1).

Storage

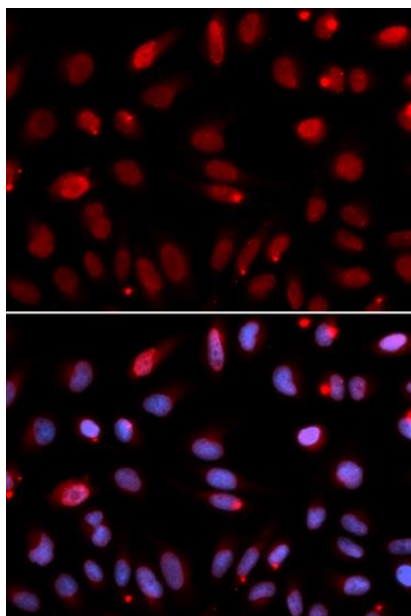
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

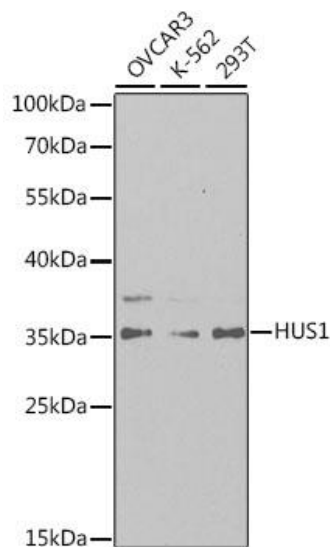
HUS1; hHUS1; checkpoint protein HUS1

Application

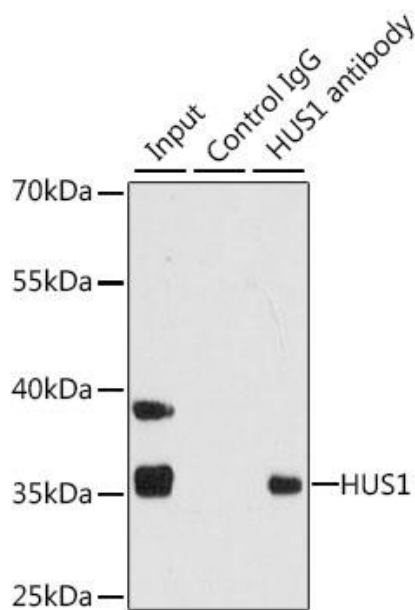
WB, IF, IP; Recommended dilution: WB 1:500 - 1:2000, IF 1:50 - 1:100, IP 1:50 - 1:200



Immunofluorescence analysis of U2OS cells using HUS1 Polyclonal Antibody. Blue: DAPI for nuclear staining.



Western blot analysis of extracts of various cell lines, using HUS1 Polyclonal Antibody at 1:1000 dilution.
Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
Lysates/proteins: 25ug per lane.
Blocking buffer: 3% nonfat dry milk in TBST.
Exposure time: 180s.



Immunoprecipitation analysis of 200ug extracts of 293T cells using 1ug HUS1 Polyclonal Antibody. Western blot was performed from the immunoprecipitate using HUS1 Polyclonal Antibody at a dilution of 1:1000.