

SETD2 Polyclonal Antibody

(Catalog # A-3720)

Background

Huntington's disease (HD), a neurodegenerative disorder characterized by loss of striatal neurons, is caused by an expansion of a polyglutamine tract in the HD protein huntingtin. This gene encodes a protein belonging to a class of huntingtin interacting proteins characterized by WW motifs. This protein is a histone methyltransferase that is specific for lysine-36 of histone H3, and methylation of this residue is associated with active chromatin. This protein also contains a novel transcriptional activation domain and has been found associated with hyperphosphorylated RNA polymerase II.

Description

SETD2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

Q9BYW2

Purification

Affinity Purified

Immunogen

Recombinant Protein of Human SETD2

Storage

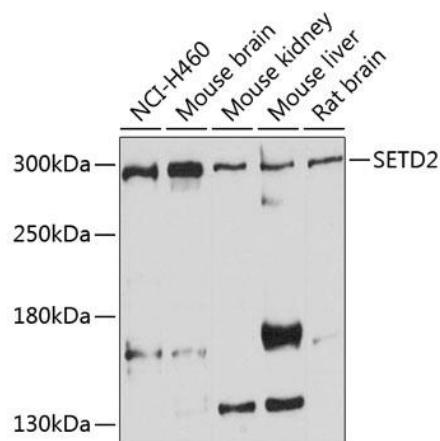
Shipped at 4°C. Store at -20°C. Avoid multiple freeze/thaw cycles.

Alternative Names

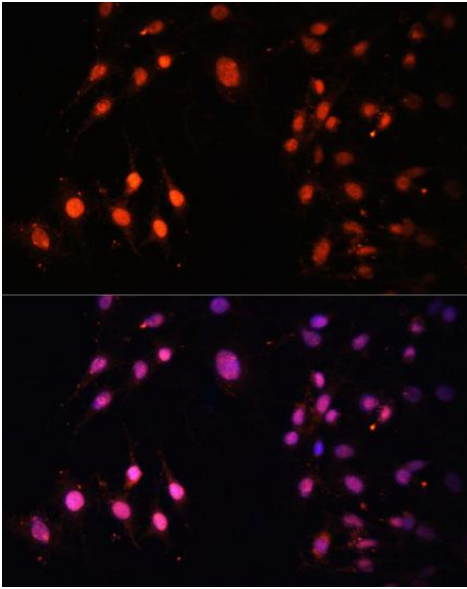
HYPB, SET2, HIF-1, HIP-1, KMT3A, HBP231, HSPC069, p231HBP

Application

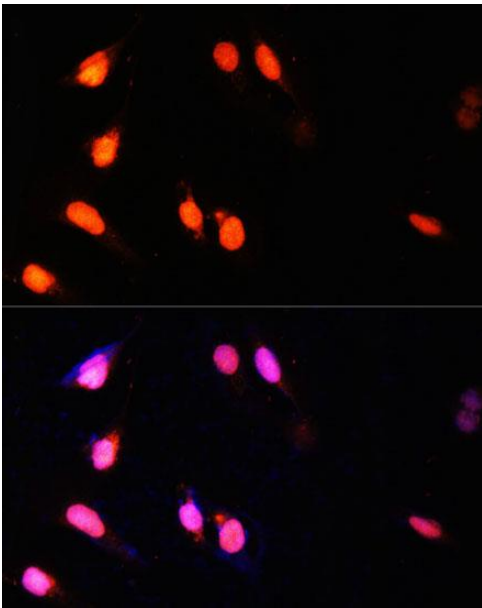
WB, IF; Recommended dilution: WB 1:500 - 1:2000; IF 1:50 - 1:200



Western blot analysis of extracts of various cell lines, using SETD2 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.



Immunofluorescence analysis of C6 cells using SETD2 Polyclonal Antibody at dilution of 1:100 (40x lens). Blue: DAPI for nuclear staining.



Immunofluorescence analysis of U-2 OS cells using SETD2 Polyclonal Antibody at dilution of 1:100 (40x lens). Blue: DAPI for nuclear staining.