
SUZ12 Polyclonal Antibody

(Catalog # A73579)

Background

This zinc finger gene has been identified at the breakpoints of a recurrent chromosomal translocation reported in endometrial stromal sarcoma. Recombination of these breakpoints results in the fusion of this gene and JAZF1. The protein encoded by this gene contains a zinc finger domain in the C terminus of the coding region.

Description

SUZ12 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

Q15022

Purification

Affinity Purified

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 470-739 of human SUZ12 (NP_056170.2).

Storage

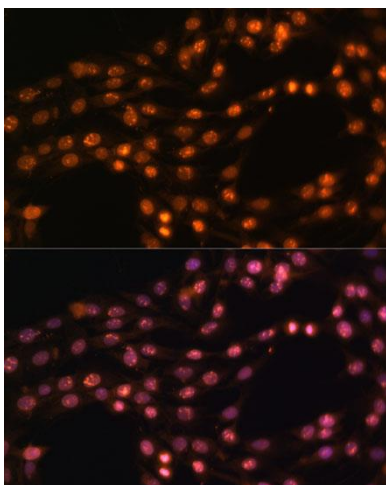
Shipped at 4°C. Store at -20°C. Avoid freeze / thaw cycles.

Alternative Names

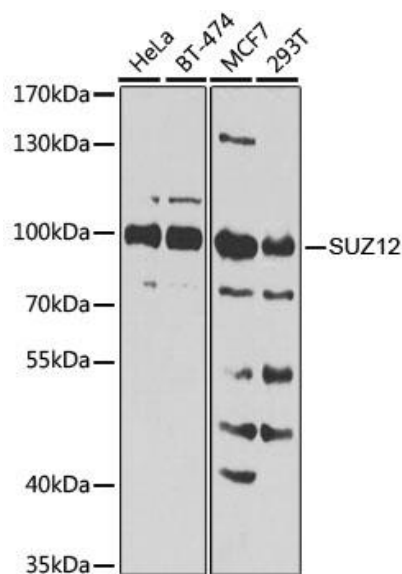
SUZ12; CHET9; JJAZ1

Application

WB, IF, IP Recommended dilution: WB 1:500 - 1:2000, IF/ICC 1:50 - 1:100, IP 1:50 - 1:100



Immunofluorescence analysis of C6 cells using SUZ12 Polyclonal Antibody at dilution of 1:100 (40x lens). Blue: DAPI for nuclear staining.

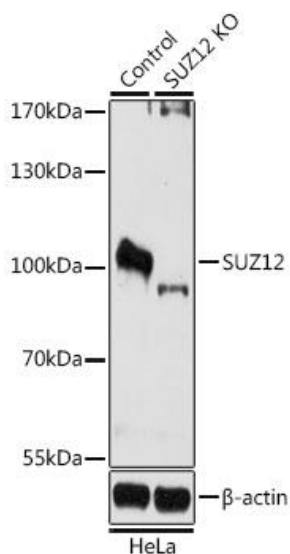


Western blot analysis of extracts of various cell lines, using SUZ12 Polyclonal Antibody at 1:1000 dilution.

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: 90s.

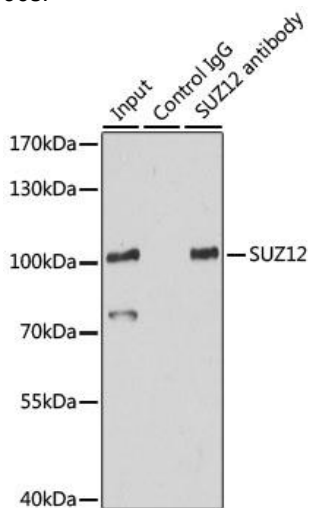


Western blot analysis of extracts from normal (control) and SUZ12 knockout (KO) HeLa cells, using SUZ12 Polyclonal Antibody at 1:1000 dilution.

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: 90s.



Immunoprecipitation analysis of 200ug extracts of HeLa cells using 3ug SUZ12 Polyclonal Antibody. Western blot was performed from the immunoprecipitate using SUZ12 Polyclonal Antibody at a dilution of 1:500.