

NSD2 / MMSET Antibody

Catalog No. 13-0002
Lot No. 15096002
Pack Size 100 µg



Type Polyclonal **Host** Rabbit
Mol. Wgt. 180 kDa **Reactivity** Human
Format Aff. Pur. IgG **Appl.** ChIP, IF, WB

Product Description:

Affinity purified rabbit polyclonal antibody recognizing human NSD2 for use in ChIP, IF and Western blot. NSD2 is a SET domain containing histone methyltransferase that catalyzes the methylation of histone H3 at lysine 36. NSD2 is also known as MMSET and WHSC1 and is implicated in several types of cancer.

Immunogen:

Recombinant protein corresponding to amino acids 959 - 1365 of human NSD2 (NM_133330).

Formulation:

200 µl immunoaffinity purified rabbit IgG at 0.5 mg/ml in PBS, pH7.5, 30% glycerol and 0.035% sodium azide.

Storage and Stability:

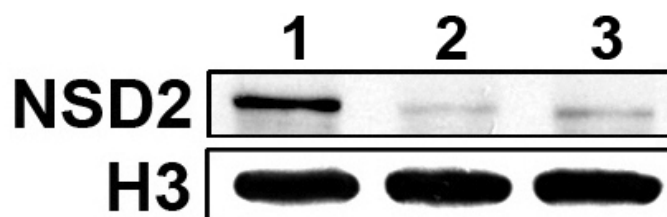
Stable for 2 years at -20°C from date of receipt.

Application Notes:

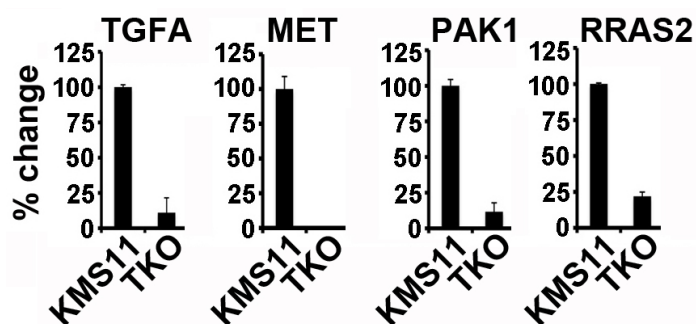
This NSD2 / MMSET antibody can be used in chromatin IP (ChIP, 2-5 µg per IP), immunofluorescence (IF, 1-2 µg/ml) and Western blotting (0.2 - 1 µg/ml).

References Using this Product:

Kuo AJ et al (2011). *Mol Cell* 44: 609-620.



Western Blot Data: Western blot using NSD2 antibody (1 µg/ml) or a total histone H3 antibody as a control. Lane 1: 100 ng KMS11 whole cell extract. Lanes 2 and 3: Extracts from KMS11 cells in which NSD2 was knocked down with RNAi.



ChIP Data: NSD2 antibody tested by ChIP (4 µg per ChIP) using chromatin from 1x10⁷ KMS11 cells or cells with the NSD2 gene knocked out (TKO). Real-time qPCR was performed on DNA purified from ChIP reactions using gene body-specific PCR primers for each gene as indicated, normalized to wild type.

Applications Key: ChIP: Chromatin IP; ChIP-seq: Chromatin IP sequencing; E: ELISA; FACS: Flow cytometry; IF: Immunofluorescence; IHC: Immunohistochemistry; IP: Immunoprecipitation; WB: Western Blotting

Reactivity Key: B: Bovine; Ce: *C. elegans*; Ch: Chicken; Dm: *Drosophila*; Eu: Eukaryote; H: Human; M: Mouse; Ma: Mammal; R: Rat; Sc: *S.cerevesiae*; Sp: *S. pombe*; WR: Wide Range (predicted); X: Xenopus; Z: Zebrafish