

Dimethyl-Arginine, asymmetric antibody (pAb)

Catalog No: 39231, 39232

Isotype: Serum

Application(s): ChIP, ChIP-Seq, ICC, IF, WB

Reactivity: Other (Not Species Specific)

Purification: None

Host: Rabbit

Volume: 200 µl, 10 µl

Background: Dimethyl-Arginine, asymmetric (or Asymmetric dimethylarginine, ADMA) interferes with the production of nitric oxide (NO). Arginine methylation is an important post-translational modification catalyzed by the PRMT/CARM family of methyltransferases.

Immunogen: This Dimethyl-Arginine, asymmetric antibody was raised against peptide containing dimethyl-arginine (asymmetric).

Buffer: Rabbit serum containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Validated Applications:

ChIP: 10 µl per ChIP

ChIP-Seq: 10 µl each

ICC/IF: 1:250 - 1:500 dilution

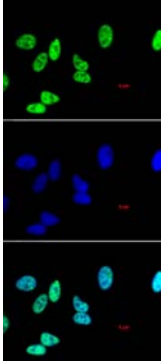
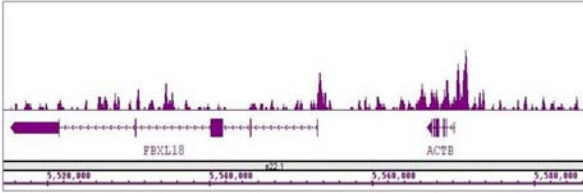
WB: 1:500 - 1:2,000 dilution

Storage and Guarantee: Antibodies in solution can be stored at -20°C for 2 years. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.

Dimethyl-Arginine, asymmetric antibody (pAb) tested by ChIP-Seq.

ChIP was performed using the ChIP-IT[®] High Sensitivity Kit (Cat. No. 53040) with 30 µg of chromatin from HeLa cells and 10 µl of antibody. ChIP DNA was sequenced on the Illumina HiSeq and 13 million sequence tags were mapped to identify Dimethyl-Arginine, asymmetric binding sites. The image shows binding across a region of chromosome 7. You can view the complete data set in the UCSC Genome Browser, starting at this specific location, here.



.Dimethyl-Arginine, asymmetric (pAb) tested by immunofluorescence.

Staining of HeLa cells with Dimethyl-Arginine, asymmetric pAb (1:250 dilution, top panel) and DAPI (middle panel), and a merge of both images (bottom panel).

Dimethyl-Arginine, asymmetric (pAb) tested by Western blot.

Detection of Dimethyl-Arginine by Western blot. The analysis was performed using 20 µg HeLa nuclear extract and Dimethyl-Arginine, asymmetric pAb at a 1:1,000 dilution

