

Asymmetric Dimethyl-arginine Antibody (ADMA)

Asym26

Catalog No. 13-0011
Lot No. 13281001
Pack Size 100 µl

Type Polyclonal **Host** Rabbit
Mol. Wgt. N/A **Reactivity** H, M, WR
Format Serum **Appl.** WB

Product Description:

Asymmetric dimethylation of arginine (ADMA) is a post-translational modification catalyzed by type I arginine methyltransferase enzymes and found on many proteins, including RNA binding proteins and histones. Asym26 recognizes ADMA present at GAR (glycine-arginine rich sequences).

Immunogen:

Synthetic peptide KFGGRGGGRGGGRGGFGGRGGRG with arg residues containing ADMA, conjugated to KLH.

Formulation:

Rabbit serum with 30% glycerol and 0.035% sodium azide.

Storage and Stability:

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

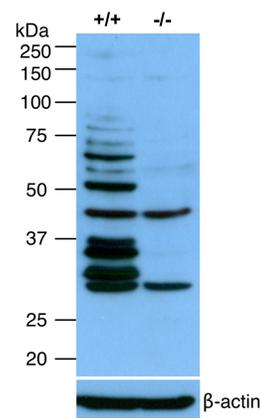
Application Notes:

Asym26 antibody is useful for Western blotting (1:500 - 1:2,000 dilution) to detect ADMA on a variety proteins.

References Using this Product:

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



Western Blot Data: Western blot using Asymmetric Dimethyl-arginine Antibody (ADMA) Asym26 antibody (1:500 dilution) on mouse embryonic fibroblast (MEF) whole cell extract derived from cells with (+/+) or without (-/-) the PRMT1 gene. Asym26 detects multiple proteins in the “PRMT1 +/+” cell extract lane, indicating proteins containing ADMA. PRMT1 is responsible for the majority of ADMA in mammals, and thus the majority of