

# Symmetric Dimethyl-arginine Antibody (SDMA)

## Sym10

**Catalog No.** 13-0012  
**Lot No.** 14125002  
**Pack Size** 100 µl

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

### Product Description:

Symmetric dimethylation of arginine (SDMA) is a post-translational modification catalyzed by type II arginine methyltransferase enzymes and found on many proteins, including those involved in transcriptional regulation, mRNA splicing, DNA repair and nuclear transport. Sym10 recognizes SDMA present at GAR (glycine-arginine rich sequences).

### Immunogen:

KLH-conjugated peptide with the sequence KRGRGRGRG in which arg residues were symmetrically dimethylated.

### 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:

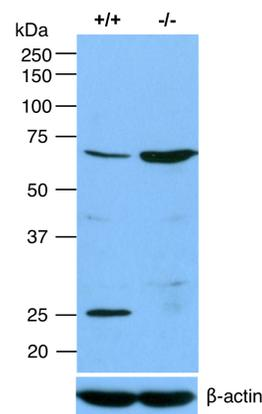
Sym10 antibody is useful for Western blotting (1:500 - 1:2,000 dilution) to detect SDMA on a variety proteins.

### References Using this Product:

Boisvert FM et al (2002). *J Cell Biol* 159: 957-969.  
Boisvert FM et al (2003). *Mol Cell Proteom* 2: 1319-1330.  
Côte J et al (2005). *J Biol Chem* 280: 28476-28483.  
Deng X et al (2010). *PNAS USA* 107: 19114-19119.  
Jung GA et al (2011). *Exp Mol Med* 43:550-560.

**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. cerevisiae*; Sp: *S. pombe*; WR: Wide Range (predicted); X: Xenopus; Z: Zebrafish



**Western Blot Data:** Western blot using Symmetric Dimethyl-arginine Antibody (SDMA) Sym10 (1:500 dilution) on mouse embryonic fibroblast (MEF) whole cell extract derived from cells with (+/+) or without (-/-) the PRMT5 gene. Sym10 detects multiple proteins in the "PRMT5 +/+" cell extract lane, indicating proteins containing SDMA. PRMT5 is responsible for the majority of SDMA in mammals. The relative migration