

Anti-Methylated Arginine (N^G-Mono-methyl)

Rabbit Polyclonal Antibody

Catalog # A95-67R

Lot # E360-2

Cited Applications

For WB(1:1000), ELISA (1:2000), IP (5ng/µg protein sample), IHC (5ng/µl) and IF (5ng/µl).

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes proteins with N^G-mono-methylated arginine residues.

Cross Reactivity

Pan-specific antibody. No reaction to acetylated proteins, di-methylated and tri-methylated proteins.

Host/Isotype/Clone#

Rabbit

Immunogen

N^G-mono-methylated arginine-KLH conjugates

Formulation

PBS, 50% glycerol, pH7.

Stability

Store at 4°C (add 0.1% NaN₃) for several months, and at -20°C for longer periods. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles.

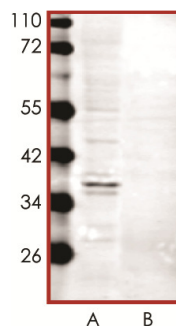
Scientific Background

Post-translational modification of proteins and peptides is a robust way to regulate function of existing proteins or peptides. Methylation on Arginine residues is one example of post-translational modification and is performed by a variety of protein methyl transferases in the cell. Arginine methylation of Rad9 for example is performed by protein arginine methyltransferase 5 (PRMT5) is important for regulation of Rad9 function and is a major element for maintaining genome integrity (1). The Methylated Arginine antibody detects Methylation on Arginine residues in proteins and peptide.

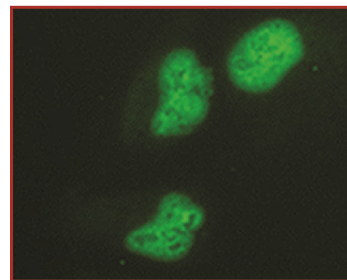
References

1. He, W. et al: A role for the arginine methylation of Rad9 in checkpoint control and cellular sensitivity to DNA damage. *Nucleic Acids Res.* 2011 Feb 14.

Sample Data



Representative western blot of the mono-methylated arginine protein profile with Anti-Methylated Arginine (N^G-Mono-methyl) (1:1000) using human melanoma (MMRU) cell lysate (A) and the MMRU lysate plus additional mono-methyl arginine peptide inhibitor (B).



Immunofluorescent stain of human melanoma (MMRU) cell with Anti-Methylated Arginine (N^G-Mono-methyl).

Anti-Methylated Arginine (N^G-Mono-methyl)

Rabbit Polyclonal Antibody

Catalog Number A95-67R

Specific Lot Number E360-2

Purification

By affinity chromatography on a mono-methyl arginine agarose column

Concentration

0.25 µg/µl

Stability

1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.