

Anti-5-Methylated Cytosine (5-methyl cytosine)

Sheep Polyclonal Antibody

Catalog # C92-67S

Lot # O2206-1

Cited Applications

WB, IP

WB: 1:500,000

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

Specificity

Recognizes the 5-Methylated Cytosine (5-methyl cytosine)

Host/Isotype/Clone#

Sheep, IgG

Immunogen

5-Methylated Cytosine (5-methyl cytosine)

Formulation

PBS

Stability

1yr at -20°C from date of shipment

Anti-5-Methylated Cytosine (5-methyl cytosine)

Rabbit Polyclonal Antibody

Catalog Number C92-67S

Specific Lot Number O2206-1

Purification Affinity chromatography

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.