



6405 Mira Mesa Blvd, Suite 100

San Diego, CA 92121

Tel: 1.858.202.1401

Fax: 1.858.481.8694

Email: support@bpsbioscience.com

Data Sheet

Anti- Adenosyl Homocysteine Antibody

Rabbit polyclonal to Adenosyl Homocysteine

Catalog #: 52030 **Lot#:** 220622

Formulated in:

PBS pH 7.2, 0.09% Sodium Azide

Purification:

Ammonium sulfate fractionation

Form : Liquid

Concentration: 5 mg/ml

Stability: Shipped at 4°C. Upon delivery aliquot and store at -20°C. Avoid freeze / thaw cycles.

Relevance:

Adenosyl Homocysteine is a competitive inhibitor of S-adenosyl-L-methionine-dependant methyl transferase reactions and therefore may play a key role in the control of methylations via regulation of the intracellular concentration of adenosylhomocysteine.

Immunogen:

Adenosyl Homocysteine conjugated to KLH.

Raised In: Rabbit

Isotype: IgG

Purification: Purified IgG prepared by affinity chromatography on Protein G

Specificity / Reactivity:

Anti-Adenosyl homosysteine antibody recognizes S-Adenosyl Homocysteine, an intermediate in the synthesis of the amino acid cysteine.

Anti-adenosyl homosysteine antibody shows the following reactivities with related compounds:

S-Adenosyl Methionine 2.5%

Adenosine 0.27%

Homocysteine <0.1%

Cystathionine <0.1%

L-Cysteine <0.1%

Glutathione <0.1%

Tested Applications: ELISA

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: support@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com