



Mouse monoclonal antibody to S-Adenosylhomocysteine (Clone 301-3) (50µl)

Catalog Number:

SAH13-A50

For Research Use Only. Not for use in diagnostic procedures.

v. 1.0

Eagle Biosciences, Inc.

20A Northwest Blvd., Suite 112, Nashua, NH 03063

Phone: 617-419-2019 Fax: 617-419-1110

www.EagleBio.com

Properties

- **Form:** Liquid
- **Storage instructions:** Store at 4°C, -20°C for long term storage
- **Storage buffer:** PBS 10mM pH 7.4 (NaCl 150mM), Sodium azide 0.02%, BSA 10mg/ml or PBS 10mM pH 7.4 (NaCl 150mM), Sodium azide 0.02%, Glycerol 50%, BSA 10mg/ml
- **Purity:** >95% Purified from mouse ascites fluid by affinity chromatography
- **Clonality:** Monoclonal
- **Clone number:** 301-3
- **Immunoglobulin isotype:** IgG3
- **Affinity:** $K_a = 8.32 \times 10^8 \text{ L/mol}$ ($1.20 \times 10^{-9} \text{ M}$)
- **Specificity:** Shows the following reactivities with related compounds:
SAdenosylhomocysteine: 100%, S-Adenosylmethionine: < 3%, Adenosine: <1%, Homocysteine: <1%, L-Cysteine: <1%, Glutathione: <1%, L-Cystathionine: <1%.
- **Immunogen:** S-Adenosylhomocysteine conjugated to BSA

Applications

The application notes include recommended starting dilutions. Optimal dilutions/concentrations should be determined by the end user. Higher dilution than suggested maybe used in IHC and IF. The product may be used in other not-yet-tested applications.

Application	Notes
cELISA	1:4000/8000
FCM	1:200
IHC	1:200

Target

S-adenosylhomocysteine is a competitive inhibitor of S-adenosylmethionine-dependant methyl transferase reactions. Therefore, it plays a key role in the control of methylation via regulation of the intracellular concentration of S-adenosylhomocysteine.

Cellular localization Cytoplasm, nuclear



Anti-Adenosylhomocysteine antibody [301-3]

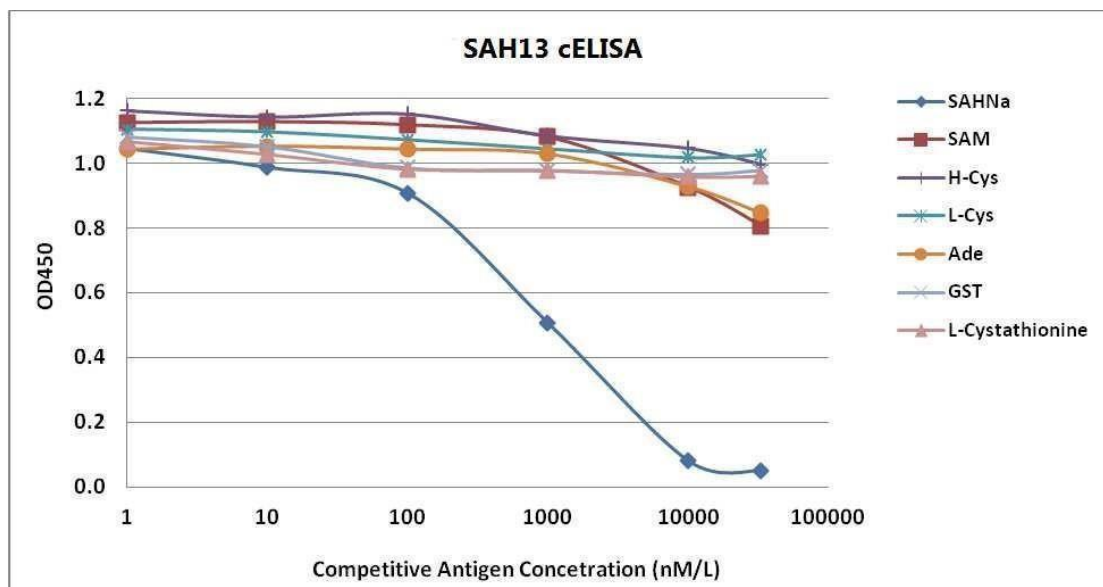


Figure 1 Competitive ELISA using anti-S-Adenosymethionine monoclonal antibody [301-3]
The 0.5 µg/ml of SAH-BSA was coated into 96 wells. Serial dilution of SAH standard (SAHNa), SAdenosylmethionine, Homocysteine (H-Cys), L-Cysteine (L-Cys), Adenosine (Ade), Glutathione (GST), L- Cystathionine (L-CTT) and properly diluted MA00303 were added. HRP conjugated Goat anti-Mouse IgG antibody was used to develop the color. OD450 value was measured on each well.

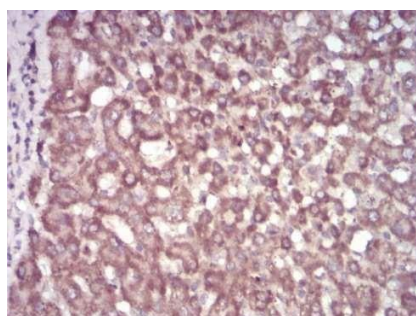


Figure 2 Immunohistochemistry staining was performed using the antibody with benign liver tissue adjacent to carcinoma. Brown areas indicated strong positive staining in nuclear and cytoplasmic areas.

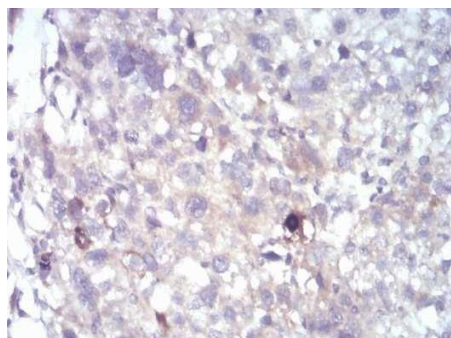


Figure 3 The same samples as in Figure 2 from liver cancer area. Cytoplasmic and nuclear areas showed negative staining.

Warranty Information

Eagle Biosciences, Inc. warrants its Product(s) to operate or perform substantially in conformance with its specifications, as set forth in the accompanying package insert. This warranty is expressly limited to the refund of the price of any defective Product or the replacement of any defective Product with new Product. This warranty applies only when the Buyer gives written notice to the Eagle Biosciences within the expiration period of the Product(s) by the Buyer. In addition, Eagle Biosciences has no obligation to replace Product(s) as result of a) Buyer negligence, fault, or misuse, b) improper use, c) improper storage and handling, d) intentional damage, or e) event of force majeure, acts of God, or accident.

Eagle Biosciences makes no warranties, either expressed or implied, except as provided herein, including without limitation thereof, warranties as to marketability, merchantability, fitness for a particular purpose or use, or non-infringement of any intellectual property rights. In no event shall the company be liable for any indirect, incidental, or consequential damages of any nature, or losses or expenses resulting from any defective product or the use of any product. Product(s) may not be resold, modified, or altered for resale without prior written approval from Eagle Biosciences, Inc.

For further information about this kit, its application or the procedures in this kit, please contact the Technical Service Team at Eagle Biosciences, Inc. at info@eaglebio.com or at 866-411-8023.