



EAGLE
BIOSCIENCES

**Horseradish peroxidase (HRP)
conjugated anti-SAH monoclonal
antibody
(Clone 301-3)**

Catalog Number:

SAHH3-A50 50 μ l

SAHH3-A100 100 μ l

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 -20°C
- **Storage buffer:** PBS 10mM pH 7.4 (NaCl 150mM), Proclin 0.1%, Glycerol 50%, BSA 10mg/ml
- **Purity:** >95% Purified from mouse ascites fluid by affinity chromatography
- **Clonality:** Monoclonal
- **Clone number:** 301-3
- **Immunoglobulin isotype:** IgG3
- **Specificity:** Shows the following reactivities with related compounds:
SAdenosylhomocysteine: 100%, S-Adenosylmethionine: ~1.5%, 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
-------------	-------

eLISA

1/2,000 – 1/6,000 (depending on the amount of coating)

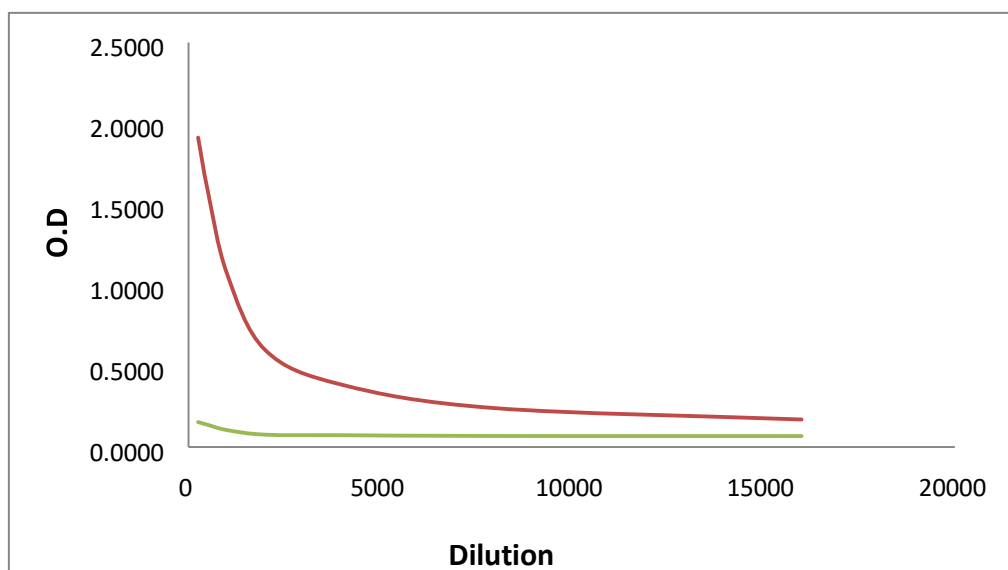




Figure 1. Titer of HRP-301-3 The 0.5 $\mu\text{g/ml}$ of SAH coating standard was coated on micro-titer wells (The top curve). The HRP-301-3 was serially diluted and added to the wells. After incubation of about 60 minutes, substrates were added and OD450 was measured. The flat curve in the bottom showed the experiment control where no SAH was coated to the plate.

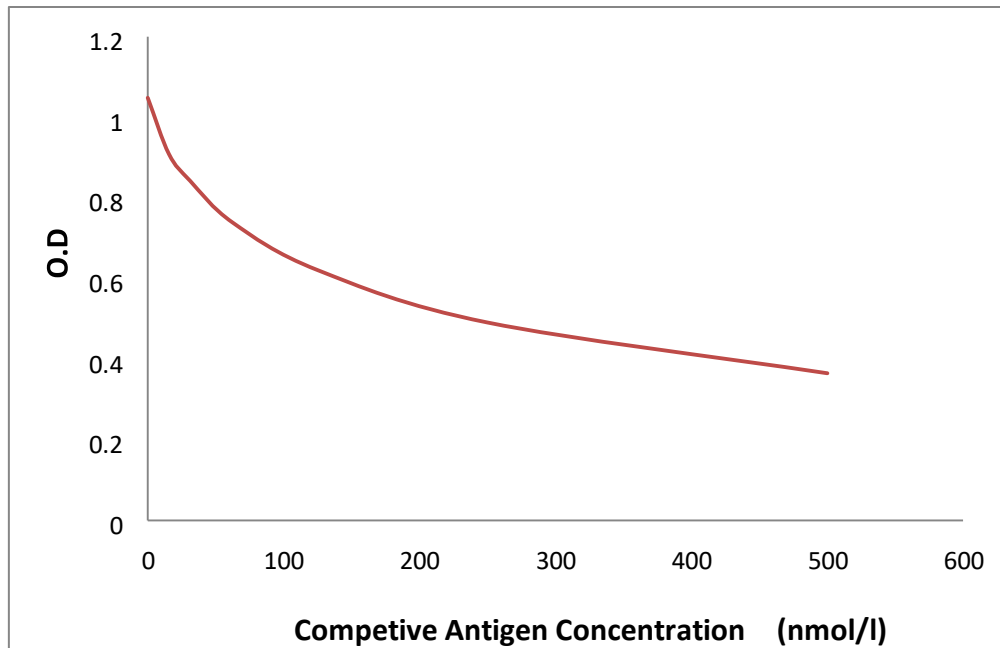


Figure 2. The standard curve with HRP-301-3 in direct competitive ELISA. The 0.5 $\mu\text{g/ml}$ of SAH coating standard was coated on micro-titer wells. Serial dilution of SAH standard was added along with 1:1,500 diluted HRP-301-3 and was incubated for about 60 minutes. The substrates were then added and OD450 was measured.

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. SAH also plays an important role in other one-carbon metabolic pathways.

Cellular localization Cytoplasm, nucleus



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.