



**EAGLE**  
BIOSCIENCES

## **Biotin Conjugated SAH (S-Adenosylhomocysteine) (6- aminohexanoic acid)**

Catalog Number:

SAH6A-A50 50µg, 25µl

SAH6A-A100 100µg, 50µl

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

*v. 1.0*

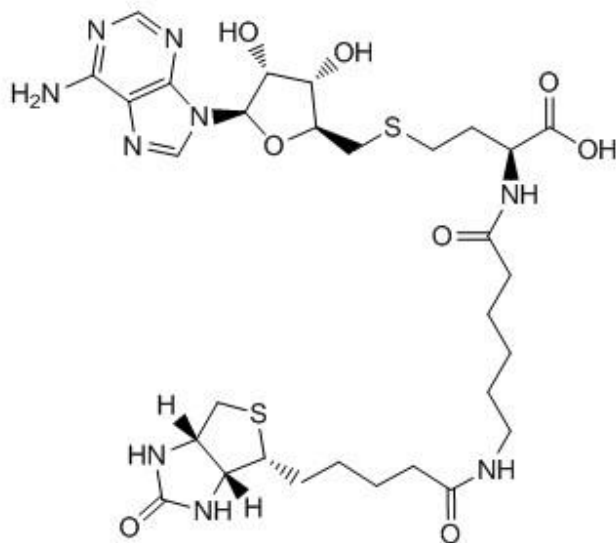
---

Eagle Biosciences, Inc.  
20A Northwest Blvd., Suite 112, Nashua, NH 03063  
Phone: 617-419-2019 Fax: 617-419-1110  
[www.EagleBio.com](http://www.EagleBio.com)



## Properties

- **Form:** Liquid
- **Molecular formula:**  $C_{30}H_{45}N_9O_8S_2$
- **Molecular Weight:** 723.86
- **Structure:**

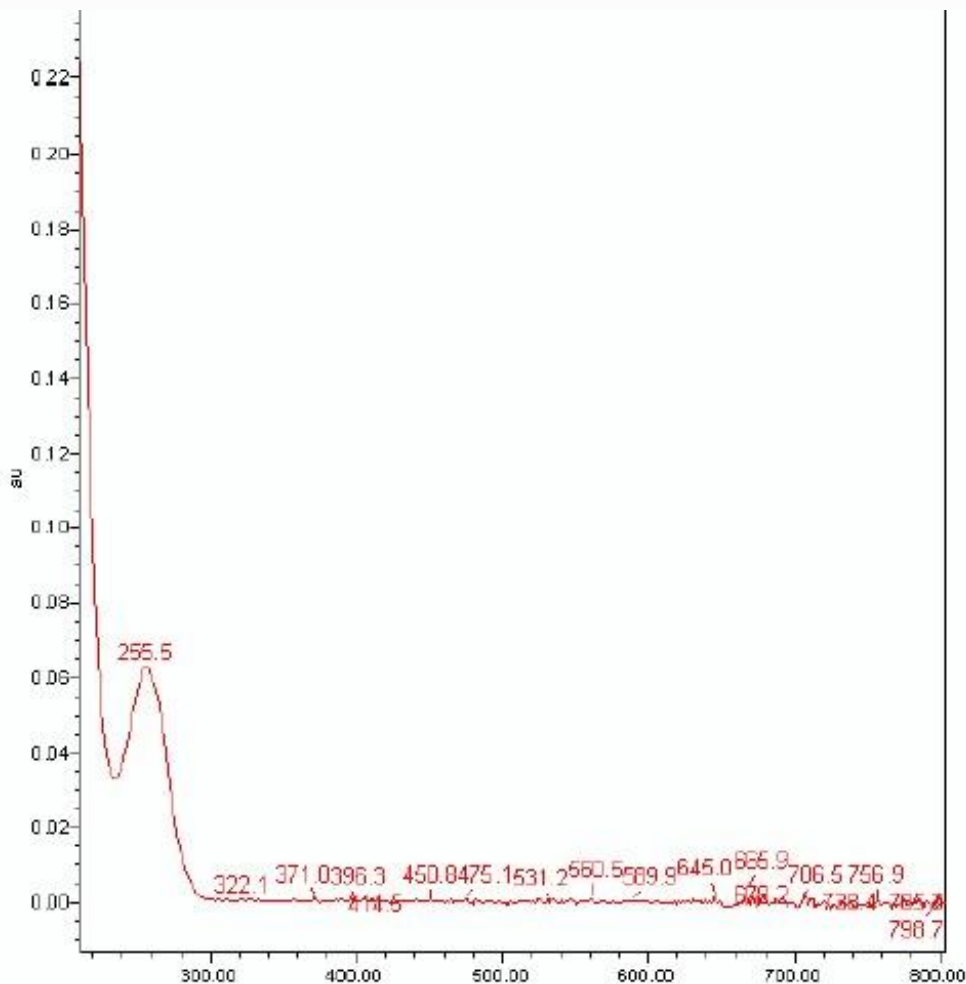


- **Storage instructions:** Store at  $-20^{\circ}\text{C}$
- **Storage buffer:** PB 20mM pH 7.4, BSA 10 mg/ml, Proclin 0.1%
- **Concentration:** 2 mg/ml
- **Purity:** 100% (HPLC)



### Verification

The ninhydrin colorimetry method was used to prove the conjugated products have the SAH component. Ultraviolet absorption spectrum of the conjugated products showed peaks at 255.5nm. SAH UV absorption spectrum does not show any peak.



### Applications

The use of this antibody in the following application has been tested. Optimal concentrations should be determined by the end user. The product may be used in other not-yet-tested applications.

Application	Notes
<b>Competitive ELIS</b>	No difference between Biotin-conjugated and unconjugated SAH molecules in their capabilities to competitively bind anti-SAH antibody.



## **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](mailto:info@eaglebio.com) or at 866-411-8023.