



Biotinylated BSA

Applicable to:

2050-0020 (20mg)

Release 2

16/02/2016

Product Details

Specificity	Avidin, Streptavidin, Neutravidin, anti-biotin antibody
Host species	Bovine
Format	Lyophilized
Buffer	Citrate buffer
Preservatives	None
Shipping	Ambient temperature
Storage	-20°C
Shelf life	2 years

Shipping conditions

The product is shipped at ambient temperature. Store the product at -20°C upon receipt.

Potential Applications

Including, but not limited to:

- ELISA
- Immunoblotting
- Immunohistochemistry

Background Information

Biotin is a small molecule with a molecular weight of 244Da. It is also known as Vitamin H, and it is crucial for a number of biological processes. Biotin-binding to avidin and streptavidin is resistant to extremes of heat, pH and proteolysis, making the capture of biotinylated molecules possible in a wide variety of environments.

Biotinylated Bovine Serum Albumin (BSA) is commonly used for measuring the binding capacity of avidin or streptavidin matrices, as a reagent for amplifying biotin-avidin detection systems, or as a positive control in biotin estimation assays.

For technical enquiries get in touch with our technical support team at:

technical.enquiries@innovabiosciences.com

For further information see our website:

www.innovabiosciences.com

Limited use license

Innova Biosciences' products are offered for research purposes alone, and are not intended for human, therapeutic or diagnostic use. The purchase of this product conveys to the buyer (whether the buyer is a not-for-profit, academic or for-profit entity) the non-transferable right to use the amount of product purchased and the components of the product for in-house research. The buyer shall not sell or otherwise transfer this product, its components, or materials prepared therefrom to any third party. The buyer shall not use this product or its components for commercial purposes. For the avoidance of doubt, 'commercial purposes' means any activity by a party for consideration and includes, without limitation, use of the product or its components (i) in the manufacturing of conjugated materials (e.g. labeled antibodies), (ii) to provide a service, information or data, (iii) for therapeutic, diagnostic or prophylactic purposes, or (iv) for repackaging/resale, whether or not such product or its components are resold for use in research.

For information on purchasing a license for commercial applications contact Innova Biosciences Ltd, Business Development Office, Babraham Hall, Babraham, Cambridge, UK, CB22 3AT. Tel +44(0)1223 496170; Fax +44(0)1223 496172.