



Streptavidin – fluorescein

Applicable to:

2014-0100 (100µg)

2014-1000 (1mg)

Release 2

16/02/2016

Product Details

Specificity	Biotin
Host species	<i>E. coli</i>
Format	Lyophilized
Buffer	Phosphate buffer
Preservatives	None
Shipping	Ambient temperature
Storage	-20°C
Shelf life	2 years

Our streptavidin conjugates are produced using our Lightning-Link® bioconjugation technology, therefore providing high quality, stable covalent conjugates for use in your application.

Background Information

Streptavidin is purified from the bacterium *Streptomyces avidinii*. It is a tetrameric protein that is composed of identical subunits, with the tetramer having a molecular weight of approximately 53kDa. Each subunit binds one biotin molecule with a very high degree of specificity and affinity. The K_d for this interaction is in the region of 10^{-15} M.

Fluorescein is a synthetic organic fluorophore, with an excitation wavelength of 494nm and a maximal emission of 520nm.

Shipping conditions

The product is shipped at ambient temperature.

Store the product at -20°C upon receipt.

Potential Applications

Including, but not limited to:

- Flow cytometry
- ELISA
- Immunoblotting
- Immunohistochemistry
- Immunocytochemistry

Related Products

Our Lightning-Link® Biotin conjugation kits allow you directly conjugate your primary antibody to biotin in under 20 minutes with only 30 seconds hands-on time:

<https://www.innovabiosciences.com/antibody-labeling-kits/biotin-streptavidin/lightning-link-biotin.html>

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.