

Streptavidin - R-Phycoerythrin (R-PE)

Applicable to: 2011-0100 (100µg), 2011-1000 (1mg)

Release 3. © EXPEDION. 11/12/2018

INTRODUCTION

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

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

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.

R-Phycoerythrin (R-PE) is a 240kDa phycobiliprotein that is isolated from red algae. R-PE has a bright red-orange fluorescence with a high quantum yield, and is brighter than other conventional fluorophores and phycobiliproteins. The absorbance spectrum of R-PE covers a range of excitation wavelengths, which provides an advantage for multi-colour immunofluorescent staining or cell-sorting. The emission maximum of R-PE is at 575nm. R-PE is used for fluorescence-based detection, often in assays which require high sensitivity.

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.expedeon.com/products/protein-antibody-conjugation/lightning-link-antibody-labeling-kits/biotin-streptavidin-labeling-kits/>

For technical enquiries get in touch with our technical support team at: technical.enquiries@expedeon.com

For further information see our website: www.expedeon.com

LIMITED USE LICENSE

Use and Use Restrictions: The Products are sold, and deliverables of any services are provided for the purposes of the buyer's internal in vitro research, development or educational use only, not for in vivo, or any therapeutic or diagnostic use, nor for resale, or for providing services or any other commercial use of any kind, including without limitation, for any transfer in any form (including as part of a kit) to a third party; for analysis or reverse engineering of the Product or for manufacturing. Products should only be used in accordance with any safety data sheets, guidance or protocols that we issue from time to time and are available for download from our Site. Protective clothing should be used at all times when handling our Products. Safety datasheets relating to all Products are available for download from the Site or upon request.

Expedeon grants no other license or rights under any intellectual property in respect of Products or services deliverables and in particular grants no license to use any Product or deliverables for any commercial purposes. Sale of Products or service deliverables by us or our authorized distributors are expressly conditional upon the customer's agreement with these restrictions, which the customer gives upon placing an order for Products or deliverables. If you wish to use any Product or deliverables for any purpose other than your own internal research as described above, you will require an additional license from Expedeon. Please contact licensing@expedeon.com.