

Protein A - HRP

Applicable to:

2000-0100 (100µg) 2000-1000 (1mg)

Release 2 16/02/2016

Product Details

Specificity	Most mammalian IgGs
Host species	E.coli and Horseradish
Format	Lyophilized
Buffer	Phosphate buffer
Preservatives	None
Shipping	Ambient temperature
Storage	-20°C
Shelf life	2 years

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

Background Information

Protein A is a 42kDa surface protein of the bacterium *Staphylococcus aureus*, which binds the Fc region of many mammalian IgGs. Its high affinity for immunoglobulins is the reason that it is often used in biochemical research.

Horseradish Peroxidase (HRP) is a metalloenzyme which has many different isoforms. It utilises hydrogen peroxide to oxidize both organic and inorganic compounds, and in the presence of the appropriate substrate (e.g. TMB, DAB, ABTS) causes a color change that can be measured spectrophotometrically.

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

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

technical.enquiries@innovabiosciences.com

For further information see our website:

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.



For information on purchasing a license for commercial applications contact Innova Biosciences Ltd, Business Development Office, Babraham Hall, Babraham, Cambridge, UK, CB22 3AT. Tel

limited use label license for Lightning-Link® products.

+44(0)1223 496170; Fax +44(0)1223 496172.