



## Protein A - Alkaline Phosphatase

Applicable to:

2001-0100 (100µg)

2001-1000 (1mg)

Release 2

16/02/2016

### Product Details

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

methyumbelliferyl phosphate (MUP) may also be used.

### 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

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.

Alkaline Phosphatase is a hydrolase enzyme. It is extracted from bovine intestine and purified using multiple precipitation, fractionation, chromatography, and filtration steps. Alkaline Phosphatase is frequently employed in immunoassays, and a commonly used colorimetric substrate for it is PNPP; the product of this reaction, p-nitrophenol, can be read at 405nm. Fluorogenic substrates such as 4-

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

[technical.enquiries@innovabiosciences.com](mailto:technical.enquiries@innovabiosciences.com)

For further information see our website:

[www.innovabiosciences.com](http://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.