

Datasheet: LNK014AP

Description:	LYNX RAPID ALKALINE PHOSPHATASE ANTIBODY CONJUGATION KIT
Name:	ALK. PHOS CONJUGATION KIT
Format:	Kit
Product Type:	Conjugation Kit
Quantity:	3 CONJUGATIONS for 10µg antibody

Product Details

Applications

This product has been reported to work in the following applications. This information is derived from testing within our laboratories, peer-reviewed publications or personal communications from the originators. Please refer to references indicated for further information. For general protocol recommendations, please visit www.bio-rad-antibodies.com/protocols.

	Yes	No	Not Determined	Suggested Dilution
Conjugation	■			

We recommend that for each conjugation the user determines the best antibody:conjugate ratio.

Product Information

LYNX Rapid Alkaline Phosphatase Antibody Conjugation Kit® enables the rapid conjugation of a pre-prepared lyophilized mixture containing Alkaline Phosphatase (AP) label to an antibody or protein. Activation of proprietary reagents within the antibody-label solution results in directional covalent bonding of AP to the antibody.

The LYNX Rapid Conjugation kit® can be used to label small quantities of antibody/protein at near neutral pH, allowing a high conjugation efficiency with 100% antibody recovery.

Reagents In The Kit

3 Vials of 10ug LYNX lyophilized AP mix
1 Vial LYNX Modifier reagent
1 Vial LYNX Quencher reagent

Preparing The Antibody

The following buffer solutions are recommended for preparing the antibody:

10-50mM amine-free buffer (e.g HEPES, MES, MOPS and phosphate*) pH range 6.5-8.5, although moderate concentrations of Tris buffer (<20mM) may be tolerated.

If possible, avoid buffers containing nucleophilic components such as primary amines and thiols (e.g. thiomersal/thimerosal) since they may react with LYNX chemicals. Azide (0.02-0.1%), EDTA and common non-buffering salts and sugars have little or no effect on conjugation efficiency.

The molar ratio of antibody: AP should be 1:1, i.e. 10ug antibody to every 10ug AP. For optimal results the antibody should be at a concentration of 1mg/ml, with a maximum volume of 10ul and a maximum antibody amount of 10ug. Antibody at a concentration of greater than 1mg/ml requires dilution. Antibody below 1mg/ml can still be used as long as the maximum volume is not exceeded. Using less than the recommended amount of antibody may result in unbound label, but this will be removed during subsequent application wash steps. Antibody below 0.5mg/ml should be concentrated before use with the kit.

* Phosphate may inhibit alkaline phosphatase activity. This kit can be used with antibodies prepared in phosphate buffer as long as the buffer will be washed away during an immunoassay procedure. Otherwise we recommend using a non-phosphate containing buffer for labeling antibodies with alkaline phosphatase.

-
- | | |
|-----------------------------|---|
| Instructions For Use | <ol style="list-style-type: none"> 1. To the antibody sample add 1ul of the Modifier reagent for every 10ul of antibody and mix gently. 2. Pipette the mixed antibody-modifier sample directly onto the LYNX lyophilized mix and gently pipette up and down twice to resuspend. 3. Replace cap onto vial and incubate at room temperature (20-25°C) for 3 hours, or overnight if preferred. 4. After incubation, add 1ul of Quencher reagent for every 10ul of antibody used. Leave to stand for 30 minutes before use. |
|-----------------------------|---|
-

- | | |
|-------------------|--|
| References | <ol style="list-style-type: none"> 1. Pérez Ruiz, E. <i>et al.</i> (2016) Evaluation of different strategies for magnetic particle functionalization with DNA aptamers. N Biotechnol. Jun 15. pii: S1871-6784(16)32260-9. [Epub ahead of print] |
|-------------------|--|
-

Storage	Store kit at -20°C only. Newly-conjugated antibody can be stored at 4°C. For long term storage however, the addition of a preservative is recommended. Storage in frost-free freezers is not recommended. This product should be stored undiluted. Avoid repeated freezing and thawing.
----------------	--

Shelf Life	12 months after despatch
-------------------	--------------------------

Health And Safety Information	Material Safety Datasheet documentation #10544 #10547 #10548 available at: Lyophilized Alkaline Phosphatase Mix (10544): https://www.bio-rad-antibodies.com/uploads/MSDS/10544.pdf Modifier Reagent (10547): https://www.bio-rad-antibodies.com/uploads/MSDS/10547.pdf Quencher Reagent (10548): https://www.bio-rad-antibodies.com/uploads/MSDS/10548.pdf
--------------------------------------	--

Licensed Use	These products and the methodology of conjugation are patent protected under United Kingdom patent number 2446088 and associated international patent applications. The purchase of this product conveys to the buyer the limited, non exclusive non-transferable right (without the right to resell repackage or further sublicense) under these patents to use the product to make conjugates for research and development purposes only. The purchaser cannot sell or otherwise transfer this product, or its components, or materials or data made using this product, or its components to a third party. Further information on purchasing licenses for diagnostic and other uses may be obtained by contacting Bio-Rad, at. Endeavour House, Langford Business Park, Langford Lane, Kidlington, Oxon. OX5 1GE UNITED KINGDOM. Tel: +44 1865 852 700. E-mail: antibodies@bio-rad.com
---------------------	---

Regulatory	For research purposes only
-------------------	----------------------------

North & South America Tel: +1 800 265 7376 Fax: +1 919 878 3751 Email: antibody_sales_us@bio-rad.com	Worldwide Tel: +44 (0)1865 852 700 Fax: +44 (0)1865 852 739 Email: antibody_sales_uk@bio-rad.com	Europe Tel: +49 (0) 89 8090 95 21 Fax: +49 (0) 89 8090 95 50 Email: antibody_sales_de@bio-rad.com
---	---	--

Printed on 11 Jul 2018

© 2018 Bio-Rad Laboratories Inc | [Legal](#) | [Imprint](#)