

Datasheet: LNK002P

Description:	LYNX RAPID HRP ANTIBODY CONJUGATION KIT
Name:	HRP CONJUGATION KIT
Format:	Kit
Product Type:	Conjugation Kit
Quantity:	3 CONJUGATIONS for 400µg 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 HRP Antibody Conjugation Kit® enables the rapid conjugation of a pre-prepared lyophilized mixture containing Horseradish peroxidase (HRP) label to an antibody or protein. Activation of proprietary reagents within the antibody-label solution results in directional covalent bonding of HRP 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 100µg LYNX lyophilized HRP 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. EDTA and common non-buffering salts and sugars have little or no effect on conjugation efficiency.

Sodium azide is an irreversible inhibitor of HRP and therefore should be avoided.

The amount of antibody used for labeling ideally should correspond to molar ratios between 1:4 and 1:1 Ab to HRP. Taking account of the molecular weights (160,000 versus 40,000), this means for that for 100µg HRP you need to add between 100-400µg of antibody. For optimal results the antibody volume should be up to 100µl, at a concentration range of 0.5-5.0mg/ml.

Instructions For Use

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

1. Bondzio, A. *et al.* (2011) Identification of differentially expressed proteins in ruminal epithelium in response to a concentrate-supplemented diet. [Am J Physiol Gastrointest Liver Physiol. 301 \(2\): G260-8.](#)
2. Lichtmanegger, J. *et al.* (2016) Methanobactin reverses acute liver failure in a rat model of Wilson disease. [J Clin Invest. 126 \(7\): 2721-35.](#)

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, although sodium azide should be avoided.
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 #10543 #10546 #10548 available at:
Lyophilized HRP Mix (10543): <https://www.bio-rad-antibodies.com/uploads/MSDS/10543.pdf>
Modifier Reagent (10546): <https://www.bio-rad-antibodies.com/uploads/MSDS/10546.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

'M317601:180705'

Printed on 11 Jul 2018