

Datasheet: LNK112PECY7

Description:	LYNX RAPID RPE-Cy7 ANTIBODY CONJUGATION KIT
Name:	RPE-Cy7 CONJUGATION KIT
Format:	Kit
Product Type:	Conjugation Kit
Quantity:	3 CONJUGATIONS for 60µg 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 RPE-Cy7 Antibody Conjugation Kit® enables the rapid conjugation of a pre-prepared lyophilized mixture containing R-Phycoerythrin (RPE)-Cy7 label to an antibody or protein. Activation of proprietary reagents within the antibody-label solution results in directional covalent bonding of RPE-Cy7 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 100ug LYNX lyophilized RPE-Cy7 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, up to 50% Glycerol and common non-buffering salts and sugars have little or no effect on conjugation efficiency.

Due to the large size of the RPE-Cy7 label, it is recommended that 50-60ug of antibody be used for every 100ug RPE-Cy7, to ensure a slight RPE-Cy7 molar excess (50ug antibody gives a 1:1 Ab:RPE-Cy7 molar ratio). For optimal results the antibody should be at a concentration of 1mg/ml, with a maximum volume of 60ul and a maximum antibody amount of 60ug. 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.

-
- | | |
|-----------------------------|--|
| Instructions For Use | <p>1.To the antibody sample add 1ul of the Modifier reagent for every 10ul of antibody and mix gently.</p> <p>2.Pipette the mixed antibody-modifier sample directly onto the LYNX lyophilized mix and gently pipette up and down twice to resuspend.</p> <p>3.Replace cap onto vial and incubate in the dark at room temperature (20-25°C) for 3 hours, or overnight if preferred.</p> <p>4.After incubation, add 1ul of Quencher reagent for every 10ul of antibody used. Leave to stand for 30 minutes before use.</p> |
|-----------------------------|--|
-

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 from date of despatch.
-------------------	----------------------------------

Acknowledgements	This product or portions thereof is manufactured under license from Carnegie Mellon University under U.S. Patent Number 5,268,486 and related patents. Cy and CyDye are trademarks of GE Healthcare Limited.
-------------------------	--

Health And Safety Information	Material Safety Datasheet documentation #10551 #10546 #10549 available at: Lyophilized RPE-Cy7 Mix (10551): https://www.bio-rad-antibodies.com/uploads/MSDS/10551.pdf Modifier Reagent (10546): https://www.bio-rad-antibodies.com/uploads/MSDS/10546.pdf Quencher Reagent (10549): https://www.bio-rad-antibodies.com/uploads/MSDS/10549.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

'M317625:180705'

Printed on 11 Jul 2018