

Datasheet: LNK203CY3

Description:	LYNX RAPID PLUS Cy3 ANTIBODY CONJUGATION KIT
Name:	Cy3 CONJUGATION KIT
Format:	Kit
Product Type:	Conjugation Kit
Quantity:	3 CONJUGATIONS for 20µg 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 Plus Cy3 antibody conjugation kit™ enables the ultra rapid conjugation of a pre-prepared lyophilized mixture containing Cyanine Dye3 (Cy3) label to an antibody or protein. Activation of proprietary reagents within the antibody-label solution results in the coupling of Cy3 to the antibody.

The LYNX Rapid Plus 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 LYNX™ Rapid Plus lyophilized Cy3 mix
1 Vial LYNX™ Rapid Plus Modifier reagent
1 Vial LYNX™ Rapid Plus 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.

Do not use buffers containing nucleophilic components e.g primary amines and thiols since they may react with LYNX™ chemicals, and Thiomersal should also be avoided. Azide (0.02-0.1%), BSA (0.1-0.5%), EDTA, Glycerol (up to 50%), and common non-buffering salts and sugars have little or no effect on conjugation efficiency.

It is recommended that 10-20ug antibody be used in each labelling reaction. For optimal results the antibody should be at a concentration of 1mg/ml, with a maximum volume of 10ul and a recommended antibody amount of 10ug. A maximum of 20ug of antibody can be used to obtain good quality conjugates as long as the maximum conjugation volume of 10ul is not exceeded. 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**
- 1.To the antibody sample add 1ul of the Rapid Plus Modifier reagent for every 10ul of antibody and mix gently.
 - 2.Pipette the mixed antibody-modifier sample directly onto the LYNX™ Rapid Plus lyophilized mix and gently pipette up and down twice to resuspend.
 - 3.Replace cap onto vial and incubate in the dark at room temperature (20-25°C) for 15 minutes, or overnight if preferred.
 - 4.After incubation, add 1ul of Rapid Plus Quencher reagent for every 10ul of antibody used, and mix gently. Leave to stand for 4 minutes before use.

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

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 #10542 #10567 #10568 available at:
 Lyophilized Cy3 Mix (10542): <https://www.bio-rad-antibodies.com/uploads/MSDS/10542.pdf>
 Rapid Plus Quencher Reagent (10567): <https://www.bio-rad-antibodies.com/uploads/MSDS/10567.pdf>
 Rapid Plus Modifier Reagent (10568): <https://www.bio-rad-antibodies.com/uploads/MSDS/10568.pdf>

Licensed Use

Bio-Rad LYNX™ conjugation kits are offered for research purposes alone, and are not intended for human, therapeutic or diagnostic use. The purchase of this conjugation kit 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. The use of this product by the buyer constitutes agreement with the terms of this limited use label license for LYNX™ products.

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

'M317641:180705'

