

## Datasheet: LNK163STR

<b>Description:</b>	LYNX RAPID STREPTAVIDIN ANTIBODY CONJUGATION KIT
<b>Name:</b>	STREPTAVIDIN CONJUGATION KIT
<b>Format:</b>	Kit
<b>Product Type:</b>	Conjugation Kit
<b>Quantity:</b>	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](http://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 Streptavidin Antibody Conjugation Kit®** enables the rapid conjugation of a pre-prepared lyophilized mixture containing Streptavidin label to an antibody or protein. Activation of proprietary reagents within the antibody-label solution results in the coupling of Streptavidin to the antibody.

LYNX Rapid Streptavidin Antibody 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 Streptavidin 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.

Do not use buffers containing nucleophilic components e.g primary amines and thiols 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.

It is recommended that 10ug 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 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.

- 
- 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.

---

**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 #10536 #10546 #10548 available at:  
Lyophilized Streptavidin Mix (10536): <https://www.bio-rad-antibodies.com/uploads/MSDS/10536.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**

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** Tel: +1 800 265 7376

**America** Fax: +1 919 878 3751

Email: [antibody\\_sales\\_us@bio-rad.com](mailto: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](mailto: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](mailto:antibody_sales_de@bio-rad.com)

'M317635:180705'

**Printed on 11 Jul 2018**