

Datasheet: STAR3000A

Description:	HISTAR DETECTION SYSTEM
Name:	HISTAR DETECTION KIT (HUMAN TISSUES ONLY)
Format:	Kit
Product Type:	Kits
Quantity:	50 TESTS

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
Immunohistology - Frozen	▪			Ready to use
Immunohistology - Paraffin	▪			Ready to use

Where this product has not been tested for use in a particular technique this does not necessarily exclude its use in such procedures. Suggested working dilutions are given as a guide only. It is recommended that the user titrates the product for use in their own system using appropriate negative/positive controls.

Intended Use

Bio-Rad HISTAR Detection System reagents are designed specifically to replace complex immunohistochemical techniques with a simple kit. This kit provides linking and labeling reagents intended for use with species specific primary antibodies for visualizing cellular antigens in tissue specimens. Visualization is accomplished using an ultra-sensitive indirect labeling method that utilizes novel polymer labeling technology. Secondary antibodies are polymerized directly with Horseradish peroxidase (HRP) into compact polymers bearing a high ratio of enzyme to antibody. This biotin-free system offers enhanced sensitivity and minimal background staining.

The HISTAR Detection System is designed for the staining of Human tissue sections. This kit is suitable for use with mouse and rabbit primary antibodies only.

All labeling and blocking reagents are pre-diluted for immediate use, and chromogen is provided in concentrated format. This product is suitable for use with the wide range of Bio-Rad ready-to-use or concentrated primary antibody reagents against human targets. Detailed instructions for use can be found in the complete Bio-Rad PDF datasheet, that can be accessed by clicking on the "Instructions For Use" link. The procedure has been briefly summarized below:

- 1) Prepare tissue sections appropriately according to the requirements of the primary antibody being used e.g. apply antigen retrieval if necessary.
- 2) Incubate paraffin-embedded sections with peroxide blocking reagent for 15 minutes at room temperature. Rinse with Phosphate Buffered Saline (PBS).
- 3) Apply serum block and incubate for 15 minutes.
- 4) Apply primary antibody and incubate for 30 minutes.
- 5) Apply reagent boost and incubate for 20 minutes. Rinse twice with PBS (5 minutes per wash).
- 6) Apply HRP polymer for 30 minutes, then rinse twice with PBS (5 minutes per wash).
- 7) Apply DAB substrate solution for 3-5 minutes, then rinse with distilled water.

[Discard unused substrate solution]

8) Counterstain and mount slides using an aqueous mounting medium such as [BUF058A](#).

Reagents In The Kit	VIAL 1, Serum Block 1 x 5ml, 5% Normal Serum in 0.01M Phosphate Buffered Saline (PBS) containing 0.1% Sodium Azide (NaN₃) VIAL 2, Reagent Boost 1 x 5ml VIAL 3, HRP Polymer 1 x 5ml VIAL 4, Peroxide Blocking Reagent 1 x 5ml, 0.03% Hydrogen Peroxide (H₂O₂) VIAL A, DAB Chromogen 1 x 2ml, 2.5% 3,3' Diaminobenzidine in stabilizing buffer VIAL B, DAB Substrate Buffer 4 x 6ml, 0.1% H₂O₂ in Citric acid buffer
Additional Reagents Required	1) Primary antibody and positive control 2) Phosphate Buffered Saline (PBS), pH7.4 (+ 0.2) 3) Counterstain 4) Mounting media Bio-Rad can supply a number of these products separately, please see the 'Recommended Useful Reagents' section.
Instructions For Use	Instructions for use can be found at www.bio-rad-antibodies.com/uploads/IFU/STAR3000A.pdf
References	1. Nishimoto, K. <i>et al.</i> (2010) Association of EMCN with Susceptibility to Rheumatoid Arthritis in a Japanese Population. J Rheumatol. 38: 221-8. 2. Dedrick, G.S. <i>et al.</i> (2011) Immunohistochemical study of human costovertebral joints: A preliminary investigation. Clin Anat. 24 (6): 741-7. 3. Alexopoulou, Z. <i>et al.</i> (2016) Deubiquitinase Usp8 regulates α-synuclein clearance and modifies its toxicity in Lewy body disease. Proc Natl Acad Sci U S A. Jul 21. pii: 201523597. [Epub ahead of print]
Storage	Store at +4°C. DO NOT FREEZE. This product should be stored undiluted.
Shelf Life	Please see label for expiry date.
Health And Safety Information	Material Safety Datasheet documentation #10233 #10362 #10282 #10456 #10185 #10090 available at: Peroxide Blocking Reagent (10233): https://www.bio-rad-antibodies.com/uploads/MSDS/10233.pdf Serum Block (10362): https://www.bio-rad-antibodies.com/uploads/MSDS/10362.pdf Reagent Boost (10282): https://www.bio-rad-antibodies.com/uploads/MSDS/10282.pdf HRP Polymer (10456): https://www.bio-rad-antibodies.com/uploads/MSDS/10456.pdf DAB Chromogen (10185): https://www.bio-rad-antibodies.com/uploads/MSDS/10185.pdf DAB Substrate Buffer (10090): https://www.bio-rad-antibodies.com/uploads/MSDS/10090.pdf
Regulatory	For research purposes only

Related Products

Recommended Useful Reagents

[10X PHOSPHATE BUFFERED SALINE \(BUF036A\)](#)

[PERMANENT AQUEOUS MOUNTING MEDIUM \(BUF058A\)](#)

North & South Tel: +1 800 265 7376

America 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

'M318898:180719'

Printed on 01 Aug 2018

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