

eBioscience™ IHC Antigen Retrieval Solution – Low pH (10X)

Catalog Number: 00-4955

For Research Use Only. Not for use in diagnostic procedures.

Product Information

Contents: eBioscience™ IHC Antigen Retrieval Solution - Low pH (10X)

REF Catalog Number: 00-4955

Formulation: 10X: 100 mM Sodium Citrate

solution, pH 6.0

Temperature Limitation: Store at room

temperature.

Batch Code: Refer to vial

LOT

Use By: Refer to vial

Description

The IHC Antigen Retrieval Solution- Low pH (10X) is designed for use during the heat induced epitope retrieval (HIER) step prior to immunohistochemistry on formalin-fixed paraffin embedded tissue sections. The use of this buffer in combination with heat (often by microwave, water bath, or pressure cooker) has been shown to restore the antigenicity of proteins modified during the formalin fixation of tissue. This buffer is supplied as a 10X stock solution. Dilution of this buffer to 1X using reagent grade water results in a 10 mM Sodium Citrate solution, pH 6.0.

Applications Reported

IHC Antigen Retrieval Solution – Low pH (10X) has been reported for use in immunohistochemical staining of formalin-fixed paraffin embedded tissue sections.

Applications Tested

IHC Antigen Retrieval Solution - Low pH (10X) has been tested by immunohistochemistry on formalin-fixed paraffin embedded tissue.

References

Lyck L. Dalmau I. Chemnitz J. Finsen B. Schrøder HD. Immunohistochemical markers for quantitative studies of neurons and glia in human neocortex. J Histochem Cytochem. 2008 Mar;56(3):201-21

Shi SR, Cote RJ, Taylor CR. Antigen retrieval techniques: current perspectives. J Histochem Cytochem. 2001 Aug;49(8):931-7.

Shi SR, Key ME, Kalra KL. Antigen retrieval in formalin-fixed, paraffin-embedded tissues: an enhancement method for immunohistochemical staining based on microwave oven heating of tissue sections. J Histochem Cytochem. 1991 Jun;39(6):741-8

Related Products

00-4956 eBioscience™ IHC Antigen Retrieval Solution – High pH (10X) 00-4960 Permount™ Mounting Medium

Not for further distribution without written consent.

Copyright © 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.