
eBioscience™ IHC/ICC Blocking Buffer - High Protein

Catalog Number: 00-4952

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

Product Information

Contents: eBioscience™ IHC/ICC Blocking Buffer - High Protein

 **Catalog Number:** 00-4952

Formulation: aqueous buffer, proteins, 0.09% sodium azide

Temperature Limitation: Store at 2-8°C.



Batch Code: Refer to vial

Use By: Refer to vial

Description

The IHC/ICC Blocking Buffer- High Protein is designed for use in immunohistochemistry and immunocytochemistry protocols which require blocking of non-specific binding sites. This blocking buffer is recommended for tissues and antibodies with high background staining, nuclear antigens, and FFPE tissue. This blocking buffer can be used both in the blocking step and as a diluent for primary and secondary antibodies during antibody incubation steps. The IHC/ICC Blocking Buffer-High Protein is compatible with organic dye-conjugated antibodies as well as unconjugated antibodies that may be used in two and three step staining protocols. This buffer is supplied as a 1X stock solution.

Applications Reported

IHC/ICC Blocking Buffer - High Protein has been reported for use in immunohistochemical staining (IHC), and immunocytochemistry (ICC).

Applications Tested

IHC/ICC Blocking Buffer - High Protein has been tested by staining with organic dye-conjugated, and unconjugated antibodies in immunohistochemistry (Frozen and FFPE) and immunocytochemistry protocols.

Related Products

00-4953 eBioscience™ IHC /ICC Blocking Buffer - Low Protein

00-4954 eBioscience™ 20X TBS Wash Buffer for IHC/ICC

00-4958 Fluoromount-G™

00-4959 Fluoromount-G™, with DAPI

44-0404 eBioscience™ StainTray™

8801-4965 eBioscience™ DAB Advanced Chromogenic Kit

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.

Tel: 888.999.1371 or 858.642.2058 • Fax: 858.642.2046 • thermofisher.com/ebioscience •

info@ebioscience.com