# invitrogen

by Thermo Fisher Scientific

# eBioscience<sup>™</sup> 20X TBS Wash Buffer for IHC/ICC

Catalog Number: 00-4954

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

## **Product Information**

Contents: eBioscience™ 20X TBS Wash Buffer for IHC/ICC

**REF** Catalog Number: 00-4954

Formulation: aqueous buffer, 0.09% sodium azide

- ł Temperature Limitation: Store at room temperature. LOT Batch Code: Refer to vial
  - Use By: Refer to vial

### Description

The 20X TBS Wash Buffer for IHC/ICC is for use in immunohistochemistry and immunocytochemistry protocols which require washing to remove non-specific binding. This wash buffer is recommended for eFluor® Nanocrystal conjugated-antibodies following antibody incubation. The TBS Wash Buffer is also compatible with organic dyeconjugated antibodies as well as unconjugated antibodies that may be used in two and three step staining protocols. The buffer is supplied as a 20X stock solution that should be diluted to 1X with distilled water.

#### **Applications Reported**

20X TBS Wash Buffer for IHC/ICC has been reported for use in immunohistochemical staining (IHC), and immunocytochemistry (ICC).

### **Applications Tested**

TBS Wash Buffer has been tested by staining with eFluor® Nanocrystal-conjugated, organic dye-conjugated, and unconjugated antibodies in immunohistochemistry (Frozen and FFPE) and immunocytochemistry protocols.

#### **Related Products**

00-4952 eBioscience™ IHC/ICC Blocking Buffer - High Protein 00-4953 eBioscience™ IHC /ICC Blocking Buffer - Low Protein 00-4958 Fluoromount-G™ 88-7071 eFluor® 605 Nanocrystal Conjugation Kit - Amine Reactive (Discontinued) 88-7072 eFluor® 650 Nanocrystal Conjugation Kit - Amine Reactive (Discontinued)

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