Assay Buffer (5X)

Catalog Number DS98200

Pub. No. MAN00018150 Rev. A.0 (30)

Product description

Volume	200 mL
Appearance/ Composition	Assay Buffer (5X): Designed to be used as (1) a blocking reagent for ELISA plates, (2) a diluent for ELISA standards and samples, (3) as a diluent for the detector antibody, and (4) as a diluent for the HRP conjugate. The Assay Buffer should appear red, slightly transparent, and non-particulate. The Assay Buffer contains a buffered solution with bovine protein as carrier and 0.5% Proclin [™] biocide as a preservative.
	Prepare a working concentration (1X) by diluting the Assay Buffer (5X) with deionized or distilled water. For example, to prepare 100 mL, add 20 mL of Assay Buffer (5x) to 80 mL of deionized or distilled water. This working Assay Buffer may be stored at 2-8°C and used for up to 2 months.
Storage and Handling	Store at 2-8°C. Please do not use the reagent beyond the expiration date printed on the label. Avoid exposure to light and extreme cold or hot temperature as this will compromise the integrity and performance of the reagent.

thermofisher.com/support | thermofisher.com/askaquestion

Manufacturer: Bender MedSystems GmbH | Campus Vienna Biocenter 2 | 1030 Vienna, Austria

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

