

5X ELISA/ELISPOT Diluent

Catalog Number 00-4202

Pub. No. MAN0018152 Rev. B.0 (31)

Product description

Description	The 5X ELISA/ELISPOT Diluent is a Phosphate Buffered Solution supplemented with bovine serum as a source of animal serum proteins. This buffer is used for blocking non-specific binding of ELISA plates, diluting recombinant standards and test samples, and making the working dilutions of the secondary detection reagent (i.e., detection antibody-biotin and AV-HRP Cat. No. 18-4100). This diluent solution also contains a preservative.
Applications Reported	Dilute the 5X ELISA/ELISPOT Diluent in deionized water reagent grade before use.
Related Products	<ul style="list-style-type: none"> • 1X TMB Solution (Cat. No. 00-4201) • Super AquaBlue ELISA Substrate (Cat. No. 00-4203)
Storage and Handling	Store at 2–8°C. 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



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

For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

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.

©2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.