invitrogen

by Thermo Fisher Scientific

eBioscience[™] Foxp3 / Transcription Factor Staining Buffer Set

Catalog Number: 00-5523

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

Product Information

REF	Contents: eBioscience [™] Foxp3 / Transcription Factor Staining Buffer Set Catalog Number: 00-5523 Handling Conditions: Use within 6 months of opening or by date indicated on the bottle		Temperature Limitation: Store at 2-8°C Batch Code: Refer to vial Use By: Refer to vial Contains sodium azide and formaldehyde
-----	--	--	--

Description

The Foxp3/Transcription Factor Staining Buffer Set has been formulated and optimized for staining with antibodies to transcription factors and nuclear proteins, such as Foxp3 and Ki-67, as well as cytokines and chemokines.

Please see the Best Protocols section for more information.

Components

Fixation/Permeabilization Concentrate (cat. 00-5123): 30 ml. Store at 2-8°C. Avoid agitation. Use within 6 months of opening. This is a 4X stock solution that must be diluted prior to use with the Fixation/Permeabilization Diluent. Dilute 1 part Concentrate with 3 parts Diluent. Caution: This solution contains formaldehyde, which is toxic and a suspected carcinogen. Contact with eyes, skin and mucous membranes should be avoided. Wear proper protective clothing and gloves.

Fixation/Permeabilization Diluent (cat. 00-5223): 100 ml. Store at 4°C. The diluent is intented to be used in combination with the Fixation/Permeabilization Concentrate.

Permeabilization Buffer (10X) (cat. 00-8333): 100 ml. Store at 2-8°C. Prior to use, this should be diluted 10-fold in distilled water. *Note: The 10X Permeabilization Buffer has a natural tendency to precipitate, however, its function is not affected by this. To clarify, the solution can be filtered after dilution to a 1X working solution.*

Applications Tested

This buffer set has been tested by intracellular staining and flow cytometric analysis of normal human peripheral blood cells or mouse splenocytes using the Best Protocols: Protocol B: One step protocol for (nuclear) intracellular proteins. This protocol may be found under the Resources Tab online.

References

Brodeur TY, Robidoux TE, Weinstein JS, Craft J, Swain SL, Marshak-Rothstein A. IL-21 Promotes Pulmonary Fibrosis through the Induction of Profibrotic CD8+ T Cells. J Immunol. 2015 Dec 1;195(11):5251-60. (**buffer**, FC, PubMed)

Berga-Bolaños R, Sharma A, Steinke FC, Pyaram K, Kim YH, Sultana DA, Fang JX, Chang CH, Xue HH, Heller NM, Sen JM.
ß-Catenin is required for the differentiation of iNKT2 and iNKT17 cells that augment IL-25-dependent lung inflammation. BMC Immunol. 2015 Oct 19;16:62. (buffer, FC, PubMed)

Related Products

12-4776 FOXP3 Monoclonal Antibody (PCH101), PE, eBioscience[™] TDS DISABLED: ABMAINT SKU (PCH101) 12-4875 EOMES Monoclonal Antibody (Dan11mag), PE, eBioscience[™] TDS DISABLED: ABMAINT SKU (Dan11mag)

12-5453 BCL6 Monoclonal Antibody (BCL-DWN), PE, eBioscience[™] TDS DISABLED: ABMAINT SKU (BCL-DWN) 12-5773 FOXP3 Monoclonal Antibody (FJK-16s), PE, eBioscience[™] TDS DISABLED: ABMAINT SKU (FJK-16s) 12-5825 T-bet Monoclonal Antibody (eBio4B10 (4B10)), PE, eBioscience[™] TDS DISABLED: ABMAINT SKU (eBio4B10 (4B10))

12-6988 ROR gamma (t) Monoclonal Antibody (AFKJS-9), PE, eBioscience™ TDS DISABLED: ABMAINT SKU (AFKJS-9)

12-9966 Gata-3 Monoclonal Antibody (TWAJ), PE, eBioscience™ TDS DISABLED: ABMAINT SKU (TWAJ) 53-5761 Nanog Monoclonal Antibody (eBioMLC-51), Alexa Fluor 488, eBioscience™ TDS DISABLED: ABMAINT

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



eBioscience[™] Foxp3 / Transcription Factor Staining Buffer Set

Catalog Number: 00-5523 For Research Use Only. Not for use in diagnostic procedures.

SKU (eBioMLC-51)

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