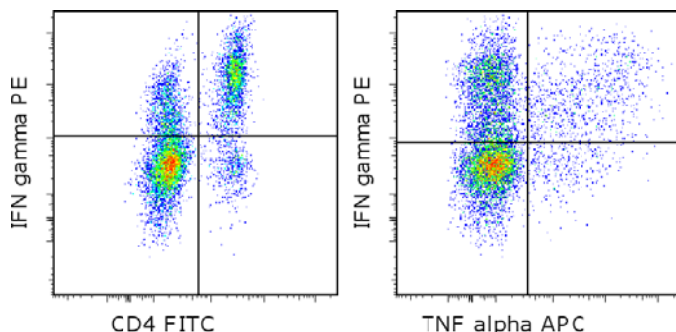


eBioscience™ Protein Transport Inhibitor Cocktail (500X)

Catalog Number: 00-4980

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



Mouse splenocytes were cultured with 1X Protein Transport Inhibitor Cocktail (500X) and 1X Cell Stimulation Cocktail (500X) (cat. 00-4970) for 16 hours. Cultures were harvested and fixed and permeabilized with IC Fixation Buffer (cat. 00-8222) and Permeabilization Buffer (cat. 00-8333). The cells were then stained with Anti-Mouse CD4 FITC (cat. 11-0042), Anti-Mouse IFN gamma PE (12-7311) and Anti-Mouse TNF alpha APC (cat. 17-7321). Cells in the lymphocyte gate were used for analysis.

Product Information

Contents: eBioscience™ Protein Transport Inhibitor Cocktail (500X)



Catalog Number: 00-4980

Concentration: 500X (Use at 1X: dilute 1/500 or to 2 uL/mL final concentration)

Handling Conditions: Wear suitable personal protective equipment and gloves.

Formulation: Brefeldin A (5.3 mM), Monensin (1 mM) in Ethanol



Temperature Limitation: Store at -20°C. Protect from light.



Batch Code: Refer to vial



Use By: Refer to vial

Description

The Protein Transport Inhibitor Cocktail (500X) is a cocktail of Brefeldin A and Monensin. It is suitable for use in assays in which prevention of protein transport to the extracellular space is desired and has been specifically formulated for use as a negative control for the Cell Stimulation Cocktail (plus protein transport inhibitors) (500X) (cat. 00-4975). Addition of this cocktail to cells in culture inhibits the intracellular protein secretory/transport pathway resulting in the accumulation of secreted proteins in the lumen of the endoplasmic reticulum and in the Golgi apparatus. This accumulation can then be detected by intracellular staining and flow cytometric analysis or western blotting.

Applications Reported

The Protein Transport Inhibitor Cocktail (500X) has been reported for use in flow cytometric analysis..

Applications Tested

The Protein Transport Inhibitor Cocktail has been tested by treatment of mouse splenocytes for 5 to 18 hours in combination with the Cell stimulation Cocktail (cat. 00-4970) followed by intracellular staining and flow cytometric analysis of IFN gamma. This cocktail is supplied as a 500X concentrate and can be directly added to culture medium at 1X (2 uL/mL).

This product is formulated with a solvent that will not freeze at the recommended storage temperatures. It is recommended that this product be kept tightly sealed and cold at all times to avoid excess evaporation.

Related Products

00-4970 eBioscience™ Cell Stimulation Cocktail (500X)
00-4975 eBioscience™ Cell Stimulation Cocktail (plus protein transport inhibitors) (500X)
00-4976 eBioscience™ Lipopolysaccharide (LPS) Solution (500X)
00-4977 eBioscience™ Phytohemagglutinin-L (PHA-L) Solution (500X)
00-4978 eBioscience™ Concanavalin A (Con A) Solution (500X)
00-8222 eBioscience™ IC Fixation Buffer

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™ Protein Transport Inhibitor Cocktail (500X)

Catalog Number: 00-4980

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

00-8333 eBioscience™ Permeabilization Buffer (10X)

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

14 March 2017 Rev. 10