# invitrogen

by Thermo Fisher Scientific

## eBioscience<sup>™</sup> Cell Stimulation Cocktail (500X)

Catalog Number: 00-4970

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



#### **Product Information**

Contents: eBioscience™ Cell Stimulation Cocktail (500X)

**REF** Catalog Number: 00-4970 Concentration: 500X (Use at 1X: dilute 1/500 or to 2 uL/mL final concentration) Handling Conditions: Wear suitable personal protective equipment and gloves. Inhibitor Cocktail (cat. 00-4980) 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

Formulation: Phorbol 12-Myristate 13-Acetate (40.5 uM), Ionomycin (670 uM) in Ethanol (500X) Temperature Limitation: Store at -20°C. Protect from light.



X

Batch Code: Refer to vial

Use By: Refer to vial

#### Description

The Cell Stimulation Cocktail (500X) is a cocktail of phorbol 12-myristate 13-acetate (PMA) and ionomycin. Treatment with PMA and ionomycin is sufficient to induce activation of many cell types to produce cytokines. This cocktail can be used to induce cytokine production in in vitro cell culture for subsequent detection in immunoassays. It can also be used in conjunction with the Protein Transport Inhibitor Cocktail (500X) (cat. 00-4980) for intracellular retention of the induced cytokine expression.

#### **Applications Reported**

The Cell Stimulation Cocktail (500X) has been reported for use in flow cytometric analysis, ELISA, and cell culture.

#### **Applications Tested**

The Cell stimulation Cocktail has been tested by treatment of mouse splenocytes for 5 to 18 hours in combination with the Protein Transport Inhibitor Cocktail (cat. 00-4980) 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 µL/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**

```
00-4975 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)
00-4980 eBioscience™ Protein Transport Inhibitor Cocktail (500X)
00-8222 eBioscience™ IC Fixation Buffer
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



### eBioscience<sup>™</sup> Cell Stimulation Cocktail (500X)

Catalog Number: 00-4970 For Research Use Only. Not for use in diagnostic procedures.

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