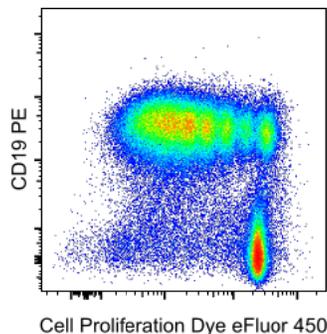


eBioscience™ Lipopolysaccharide (LPS) Solution (500X)

Catalog Number: 00-4976

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



Mouse splenocytes were labeled with 10 uM Cell Proliferation Dye eFluor® 450 (cat. 65-0842) and cultured for 3 days with Lipopolysaccharide (LPS) Solution (500X) at 2 uL per mL of culture medium. Cells were stained with Anti-Mouse CD19 PE (cat. 12-0193) and Fixable Viability Dye eFluor® 780 (cat. 65-0865). Total singlet-gated, viable cells were used for analysis.

Product Information

Contents: eBioscience™ Lipopolysaccharide (LPS) Solution (500X)

REF **Catalog Number:** 00-4976

Concentration: 500X (2.5 mg/mL)

Handling Conditions: Use in sterile environment.

Source: *Escherichia coli* 026:B6

Formulation: Sterile aqueous buffer, no sodium azide

Temperature Limitation: Store at -20°C.

LOT **Batch Code:** Refer to vial

Use By: Refer to vial

Description

The Lipopolysaccharide (LPS) Solution (500X) is a ready to use solution of LPS from *Escherichia coli* 026:B6 in aqueous solution. LPS is a major component of the cell wall of gram negative bacteria. It is highly immunogenic and strongly activates immune cells bearing the CD14/TLR4/MD2 receptor complex. This reagent is intended for use in *in vitro* culture of B cells, monocytes, macrophages and other cells that are responsive to LPS.

Applications Reported

Lipopolysaccharide (LPS) Solution (500X) has been reported for use in *in vitro* cultures.

Applications Tested

The activity of the Lipopolysaccharide (LPS) Solution (500X) has been tested by proliferation of mouse splenocytes, as measured by dilution of Cell Proliferation Dye eFluor® 450. This is a pre-titrated 500X solution and can be diluted to 2 uL per mL of culture medium. This reagent may be further-titrated for optimal performance in the assay of interest.

Under the testing conditions listed above, no change in performance is observed after 20 freeze-thaw cycles. For optimal performance, smaller aliquots may be prepared to minimize the number of freeze-thaw cycles.

References

Rittig MG, Kaufmann A, Robins A, Shaw B, Sprenger H, Gemsa D, Foulongne V, Rouot B, Dornand J. Smooth and rough lipopolysaccharide phenotypes of *Brucella* induce different intracellular trafficking and cytokine/chemokine release in human monocytes. *J Leukoc Biol.* 2003 Dec;74(6):1045-55.

Rietschel ET, Kirikae T, Schade FU, Mamat U, Schmidt G, Loppnow H, Ulmer AJ, Zähringer U, Seydel U, Di Padova F, et al. Bacterial endotoxin: molecular relationships of structure to activity and function. *FASEB J.* 1994 Feb;8(2):217-25.

Smith CI, Hammarström L, Bird AG, Kunori T, Gustafsson B, Holme T. Lipopolysaccharide and lipid A-induced human blood lymphocyte activation as detected by a protein A plaque assay. *Eur J Immunol.* 1979 Aug;9(8):619-25.

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™ Lipopolysaccharide (LPS) Solution (500X)

Catalog Number: 00-4976

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

Skidmore BJ, Chiller JM, Morrison DC, Weigle WO. Immunologic properties of bacterial lipopolysaccharide (LPS): correlation between the mitogenic, adjuvant, and immunogenic activities. J Immunol. 1975 Feb;114(2 pt 2):770-5.

Andersson J, Möller G, Sjöberg O. Selective induction of DNA synthesis in T and B lymphocytes. Cell Immunol. 1972 Aug;4(4):381-93.

Related Products

00-4505 eBioscience™ Monensin Solution (1000X)

00-4506 eBioscience™ Brefeldin A Solution (1000X)

00-4970 eBioscience™ Cell Stimulation Cocktail (500X)

00-4975 eBioscience™ Cell Stimulation Cocktail (plus protein transport inhibitors) (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)

65-0842 eBioscience™ Cell Proliferation Dye eFluor™ 450

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