

# LPS-EB

Standard lipopolysaccharide from *E. coli* O111:B4 strain; TLR4 and TLR2 ligand

Catalog code: tlrl-ebmps

<https://www.invivogen.com/lps-eb>

For research use only

Version 23F20-MM

## PRODUCT INFORMATION

### Contents

- 5 mg LPS-EB (lipopolysaccharide from *E. coli* O111:B4)
- 1.5 ml endotoxin-free water

### Storage and stability

- LPS-EB is shipped at room temperature. Upon receipt, store product at -20°C.
- Resuspended LPS-EB may be stored for 1 month at 4°C or for 6 months when aliquoted and stored at -20°C. Avoid repeated freeze-thaw cycles.

### Quality control

- Activation of TLR4 has been confirmed using HEK-Blue™ TLR4 cells.
- The presence of other bacterial components (e.g. lipoproteins) has been assessed using HEK-Blue™ TLR2 cells.

## DESCRIPTION

Lipopolysaccharide (LPS) is the principal component of Gram-negative bacteria that activates the innate immune system through its recognition by Toll-like receptor 4 (TLR4). This leads to a signaling cascade that ultimately results in the activation of NF-κB and the production of proinflammatory cytokines<sup>1</sup>. LPS-EB is a preparation of smooth (s)-form LPS purified from the Gram-negative *E. coli* O111:B4, a pathogenic serotype of *E. coli* known to cause significant gastric disease<sup>2,3</sup>.

LPS-EB is a standard lipopolysaccharide (LPS) preparation extracted by a phenol-water mixture. LPS-EB contains other bacterial components, such as lipoproteins, and therefore stimulates both TLR4 and TLR2.

1. Kuzmich N.N. *et al.*, 2017. TLR4 Signaling pathway modulators as potential therapeutics in inflammation and sepsis. *Vaccines (Basel)* 5(4):34. 2. Coleman W.G., Jr. *et al.*, 1977. Genetic analysis of *Escherichia coli* O111:B4, a strain of medical and biochemical interest. *J Bacteriol* 130:656-60. 3. Viljanen M.K. *et al.*, 1990. Outbreak of diarrhea due to *Escherichia coli* O111:B4 in schoolchildren and adults: association of Vi antigen-like reactivity. *Lancet* 336:831-4.

## PRODUCT PROPERTIES

**Species:** *Escherichia coli*

**Specificity:** TLR4 and TLR2 agonist

**Working concentration:** 10 ng-10 µg/ml

**Solubility:** 5 mg/ml in water

## METHODS

### Preparation of stock solution (5 mg/ml)

1. Add 1 ml of endotoxin-free water (provided) and homogenize.

#### Notes:

- This product can be reconstituted by injecting water through the rubber cap using a needle and syringe.
- LPS-EB stock solution may appear cloudy.

### TLR4 activation using LPS-EB

LPS-EB can be used to activate TLR4 in HEK-Blue™ TLR4 cells, that were designed to study TLR4 stimulation by monitoring NF-κB activation. Stimulation of HEK-Blue™ TLR4 cells with a TLR4 agonist activates NF-κB which induces the production of SEAP (secreted embryonic alkaline phosphatase). Levels of SEAP can be easily determined using HEK-Blue™ Detection, a cell culture medium that allows the detection of SEAP as the reporter protein is secreted by the cells.

For more information visit: <https://www.invivogen.com/hek-blue-htlr4>.

1. Add 20 µl of LPS-EB at 10 ng-10 µg/ml in a well of a 96-well plate.
2. Prepare a cell suspension ~140,000 cells per ml in HEK-Blue™ Detection.
3. Add 180 µl of the cell suspension (~25,000 cells) to each LPS-EB-containing well.
4. Incubate the plate for 6-24 h at 37°C, 5% CO<sub>2</sub>.
5. Determine SEAP levels using a spectrophotometer at 620-655 nm.

## RELATED PRODUCTS

Product	Description	Cat. Code
HEK-Blue™ Detection	SEAP Detection reagent	hb-det2
HEK-Blue™ hTLR4 Cells	Human TLR4 reporter cells	hkb-htlr4
HEK-Blue™ mTLR4 Cells	Murine TLR4 reporter cells	hkb-mtlr4
LPS-EB Ultrapure	LPS from <i>E. coli</i> O111:B4	tlrl-3pelps
LPS-SM Ultrapure	LPS from <i>S. minnesota</i>	tlrl-smmps
MPLAs	Synthetic MPLA	tlrl-mpis

## TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Asia: +852 3622-3480

E-mail: [info@invivogen.com](mailto:info@invivogen.com)