

LPS-RS Ultrapure

Ultrapure lipopolysaccharide from *Rhodobacter sphaeroides*- TLR4 antagonist

Catalog # tlr1-prslps

For research use only

Version # 15L17-MM

PRODUCT INFORMATION

Content:

- 1 mg LPS-RS Ultrapure
- 1.5 ml endotoxin-free water

Storage:

- LPS-RS Ultrapure is shipped at room temperature and should be stored at -20°C. Lyophilized product is stable 1 year at -20°C when properly stored.
- Upon resuspension, prepare aliquots of LPS-RS Ultrapure and store at 4°C for short term storage or -20°C for long term storage. Resuspended product is stable 1 month at 4°C and 6 months at -20°C. Avoid repeated freeze-thaw cycles.

BACKGROUND

Bacterial lipopolysaccharide (LPS) is the major structural component of the outer wall of all Gram-negative bacteria and a potent activator of the immune system. LPS is recognized by Toll-like receptor 4 (TLR4) which interacts with three different extracellular proteins: LPS binding protein (LBP), CD14 and myeloid differentiation protein 2 (MD-2), to induce a signaling cascade leading to the activation of NF-κB and the production of proinflammatory cytokines.

LPS consists of a polysaccharide region that is anchored in the outer bacterial membrane by a specific carbohydrate lipid moiety termed lipid A. Lipid A, also known as endotoxin, is responsible for the immunostimulatory activity of LPS. Lipid A is a glucosamine disaccharide linked to hydroxy fatty acids that are further substituted by nonhydroxylated fatty acids. The number of fatty acids is a major determinant of the immunogenicity of endotoxin. The most active form of lipid A contains six fatty acyl groups and is found in pathogenic bacteria such as *Escherichia coli* and *Salmonella* species. Underacylated lipid A structures, containing four or five fatty acids, induce markedly less host defense responses and can inhibit in a dose-dependent manner the strong endotoxic response triggered by hexa-acylated LPS.

LPS from the photosynthetic bacterium *Rhodobacter sphaeroides* (LPS-RS) is a potent antagonist of toxic LPS in both human and murine cells and also prevents LPS-induced shock in mice¹. LPS-RS is penta-acylated that appears to utilize at least two distinct mechanisms to block LPS-dependent activation of TLR4. The main mechanism consists of direct competition between under-acylated LPS and hexa-acylated LPS for the same binding site on MD-2, while the secondary mechanism involves the ability of under-acylated LPS:MD-2 complexes to inhibit hexa-acylated endotoxin:MD-2 complexes function at TLR4^{2,5}.

DESCRIPTION

LPS-RS Ultrapure is a purified preparation of the TLR4 antagonist LPS-RS (cat. code tlr1-rslps). Standard LPS-RS contains lipopeptide contaminants, and therefore stimulates Toll-like receptor 2 (TLR2), whereas LPS-RS Ultrapure only has extremely weak TLR2 activity. In order to remove TLR2 activating bacterial components, LPS-RS Ultrapure was purified by successive phenol extractions. LPS-RS Ultrapure displays strong TLR4 antagonist activity. Complete competitive inhibition of LPS activity is possible at a 100 fold excess of the antagonist. LPS-RS Ultrapure does not induce TLR4 signaling but is detected by the LAL assay, the standard endotoxin detection assay.

APPLICATION

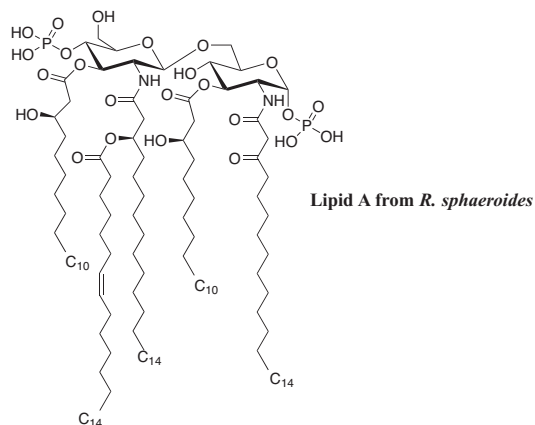
TLR4-specific inhibition

PROPERTIES

Working concentration: 10 ng - 10 µg/ml

Solubility: 1 mg/ml in water

Structure:



1. Qureshi N. *et al.*, 1999. Nontoxic RsDPLA as a potent antagonist of toxic lipopolysaccharide. H. Brade, and S. M. Opal, and S. N. Vogel, and D. C. Morrison, eds. *Endotoxin in Health and Disease* 687. Marcel Dekker, New York.
2. Coats SR. *et al.*, 2005. MD-2 mediates the ability of tetra-acylated and penta-acylated lipopolysaccharides to antagonize *Escherichia coli* lipopolysaccharide at the TLR4 signaling complex. *J Immunol.*;175(7):4490-8.
3. Teghanemt A. *et al.*, 2005. Molecular basis of reduced potency of underacylated endotoxins. *J Immunol.* 175(7):4669-76.
4. Visintin A. *et al.*, 2005. Pharmacological inhibition of endotoxin responses is achieved by targeting the TLR4 coreceptor, MD-2. *J Immunol.* 175(10):6465-72.
5. Saitoh S. *et al.*, 2004. Lipid A antagonist, lipid IVa, is distinct from lipid A in interaction with Toll-like receptor 4 (TLR4)-MD-2 and ligand-induced TLR4 oligomerization. *Int Immunol.* 16(7):961-9.

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 Hong Kong: +852 3-622-34-80

E-mail: info@invivogen.com

METHODS

Preparation of a stock solution and dilutions

- Add 1 ml endotoxin-free water (provided) to the 1 mg of LPS-RS Ultrapure and homogenize to obtain a stock solution at 1 mg/ml.
- Prepare serial dilutions using endotoxin-free water, typically from 10 ng to 10 µg/ml.

Inhibition of TLR4 activation

InvivoGen has developed a simple and convenient cell-based method to study TLR4 activation. This method relies on the use of HEK-Blue™-hTLR4 cells that coexpress human TLR4, MD2 and CD14 genes, and an optimized SEAP reporter gene placed under the control of an NF-κB-inducible promoter. Upon TLR4 activation, the cells secrete SEAP which can be readily determined using QUANTI-Blue™, a detection medium that turns blue in the presence of SEAP.

The protocol described below is performed using HEK-Blue™-hTLR4 cells, QUANTI-Blue™, the LPS from *E. coli* K12 (LPS-EK) and the antagonist LPS from *R. sphaeroides* (LPS-RS Ultrapure). Alternatively, the LPS from *E. coli* 0111:B4 (LPS-EB) can be used to activate TLR4.

Day 1:

- Prepare a HEK-Blue™-hTLR4 cell suspension at 160,000 cells/ml in DMEM supplemented with 10% heat-inactivated fetal bovine serum.

Note: Some fetal bovine serum (FBS) may contain alkaline phosphatases that can interfere with SEAP quantification. We recommend to use heat-inactivated FBS to inactivate these enzymes which are thermosensitive. Heat-inactivated FBS is available commercially or can be prepared by incubating 30 min at 56°C.

- Add 20 µl of LPS-RS Ultrapure and 160 µl of HEK-Blue™-hTLR4 cell suspension (~25,000 cells) per well of a flat-bottom 96-well plate.
- Add 20 µl of LPS-EK per well of a flat-bottom 96-well plate at the final concentrations indicated in the table below:

LPS-EK Conc. (ng/ml)	LPS-RS Conc. (ng/ml)					
	0	10	30	100	300	1000
0	0	10	30	100	300	1000
0.25	0	10	30	100	300	1000
0.5	0	10	30	100	300	1000
1	0	10	30	100	300	1000

- Incubate the plate for 18-24 h at 37°C, 5% CO₂.

Day 2:

- Prepare QUANTI-Blue™ following the instructions on the pouch.
- Add 180 µl of resuspended QUANTI-Blue™ per well of a flat-bottom 96-well plate.
- Add 20 µl of induced HEK-Blue™-hTLR4 Cells supernatant.
- Incubate the plate at 37°C incubator for 1 to 3 h.
- Determine SEAP levels using a spectrophotometer at 620-655 nm.

RELATED PRODUCTS

Product	Catalog Code
HEK-Blue™-hTLR4 Cells	hkb-htlr4
LPS-EB	tlrl-ebllps
LPS-EK	tlrl-ekllps
QUANTI-Blue™	rep-qbl

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 Hong Kong: +852 3-622-34-80
E-mail: info@invivogen.com