

LPS-RS

Lipopolysaccharide from *Rhodobacter sphaeroides*- TLR4 antagonist

Catalog # tlr1-rslps

For research use only

Version # 13106-MM

PRODUCT INFORMATION

Content:

- 5 mg *Rhodobacter sphaeroides* LPS (LPS-RS)
- 1.5 ml endotoxin-free water

Storage:

- LPS-RS 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 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.

APPLICATION

TLR4-specific inhibition in TLR2 non-responsive cells.

DESCRIPTION

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.

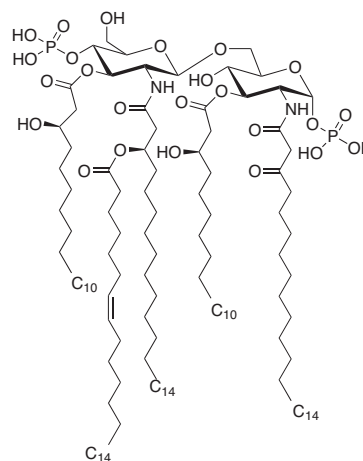
LPS consists of a polysaccharide region that is anchored in the outer bacterial membrane by a specific carbohydrate lipid moiety termed lipid A. Lipid A, also known as endotoxin, is responsible for the immunostimulatory activity of LPS. Lipid A is a glucosamine disaccharide linked to hydroxy fatty acids 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 TLR4²⁻⁵. Complete competitive inhibition of LPS activity is possible at a 100 fold excess of the antagonist. LPS-RS does not induce TLR4 signaling but is detected by the LAL assay, the standard endotoxin detection assay.

LPS-RS preparations (cat. code tlr1-rslps) contain lipopeptide contaminants, and therefore stimulate Toll-like receptor 2 (TLR2).

1. Qureshi, N., B. W. Jarvis, K. Takayama. 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.



Lipid A from *R. sphaeroides*

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 3622-3480
E-mail: info@invivogen.com

METHODS

Preparation of a stock solution and dilutions

- Add 1 ml endotoxin-free water (provided) to the 5 mg of LPS-RS and homogenize to obtain a stock solution at 5 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 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.

Note: Be aware that if your cells express TLR2, inhibition of TLR4 with LPS-RS will not be detectable, as LPS-RS activates TLR2.

The protocol described below is performed using HEK-Blue™-hTLR4 Cells QUANTI-Blue™, the ultrapure LPS from *E. coli* 0111:B4 (LPS-EB ultrapure) and the antagonist LPS from *R. sphaeroides* (LPS-RS). Alternatively, the ultrapure LPS from *E. coli* K12 (LPS-EK ultrapure) 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 FBS.

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.

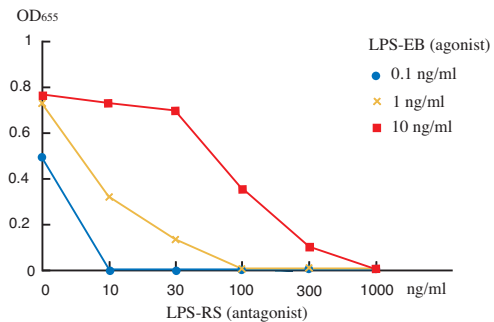
- Add 20 µl of LPS-EB and 20 µl of LPS-RS per well of a flat-bottom 96-well plate at the final concentrations indicated in the table below:

LPS-EB Conc. (ng/ml)	LPS-RS Conc. (ng/ml)					
0	0	10	30	100	300	1000
0.1	0	10	30	100	300	1000
1	0	10	30	100	300	1000
10	0	10	30	100	300	1000

- Add 160 µl of HEK-Blue™-hTLR4 cell suspension (~25,000 cells) per well.
- Incubate the plate at 37°C in a 5% CO₂ incubator for 18-24 h.

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.



Antagonistic effect of LPS-RS on LPS-stimulated TLR4: HEK-Blue™-4 cells were incubated with increasing concentrations of LPS-EB ultrapure (*E. coli* 0111:B4) and LPS-RS. After 24h incubation, TLR4 inhibition was assessed by measuring the levels of SEAP secreted in the supernatant by using QUANTI-Blue™, which turns blue in the presence of SEAP.

RELATED PRODUCTS

Product	Catalog Code
HEK-Blue™-hTLR4 Cells	hkb-htlr4
LPS-EB Ultrapure	tlr1-3pelps
LPS-EK Ultrapure	tlr1-peklps
QUANTI-Blue™	rep-qb1

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 3622-3480
 E-mail: info@invivogen.com

