

LTA-SA Purified

Purified lipoteichoic acid from *Staphylococcus aureus* - TLR2 ligand

Catalog # tlr1-pslta

For research use only

Version # 15E011-MM

PRODUCT INFORMATION

Content:

- 5 mg LTA-SA Purified
- 10 ml endotoxin-free water

Storage:

- LTA-SA Purified is shipped at room temperature and should be stored at -20 °C.
- Upon resuspension, store LTA-SA Purified at 4 °C for short-term storage or -20 °C for long storage. Avoid repeated freeze-thaw cycles. Resuspended product is stable for 1 month at 4 °C and for 6 months at -20 °C when properly stored.

Quality Control:

- The TLR2 activity has been tested using HEK-Blue™ TLR2 cells.
- The absence of other bacterial components (e.g. endotoxins) has been confirmed using HEK-Blue™ TLR4 cells.

DESCRIPTION

Lipoteichoic acid (LTA) is a major immunostimulatory component of Gram-positive bacteria. LTA is responsible for causing gram-positive sepsis. LTA shares many of the biochemical and physiological properties of the immunostimulatory lipopolysaccharide (LPS) of gram-negative bacteria. Like LPS, LTA is an amphiphile formed by a hydrophilic polyphosphate polymer linked to a neutral glycolipid. LTA stimulates immune cells through TLR2 to produce TNF- α and other inflammatory cytokines¹.

This preparation of LTA from *S. aureus* is purified following the method described by Morath *et al.*². It contains 10 times less endotoxin according to the gel clot LAL Assay than the standard preparation. At concentrations ranging from 1 ng to 10 μ g/ml, it highly activates TLR2 and no other TLRs including TLR4.

1. Schwandner R. *et al.*, 1999. Peptidoglycan- and lipoteichoic acid-induced cell activation is mediated by toll-like receptor 2. *J Biol Chem*, 274(25):17406-9. 2. Morath S. *et al.*, 2001. Structure-function relationship of cytokine induction by lipoteichoic acid from *Staphylococcus aureus*. *J.Exp.Med.* 193:393-397.

METHODS

Preparation of stock solution (5 mg/ml)

Stimulation of TLR2 can be achieved with 1 ng - 1 μ g/ml LTA-SA Purified.

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

TLR2 stimulation with LTA-SA Purified

LTA-SA Purified can be used to activate TLR2 in cells expressing this receptor such as HEK-Blue™ TLR2 cells. These cells were designed to study TLR2 stimulation by monitoring NF- κ B activation. Stimulation of HEK-Blue™ TLR2 cells with a TLR2 agonist activates NF- κ B 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: www.invivogen.com/hek-blue-trl2

- Dispense 20 μ l of LTA-SA Purified at various concentrations (1 ng - 1 μ g/ml) per well of a 96-well plate.
- Prepare a cell suspension ~280,000 cells per ml in HEK-Blue™ Detection medium and immediately add 180 μ l of the cell suspension (~50,000 cells) to each LTA-SA Purified-containing well.
- Incubate the plate for 6 - 24 h at 37 °C, 5% CO₂.
- Determine SEAP levels using a spectrophotometer at 620 - 655 nm.

RELATED PRODUCTS

Product	Catalog Code
HEK-Blue™ Detection	hb-det2
HEK-Blue™ hTLR2 cells	hkb-htlr2
Other TLR2 agonists	
FSL-1 (Synthetic diacylated lipoprotein)	tlr1-fsl
HKLM (Heat Killed <i>L. monocytogenes</i>)	tlr1-hklm
LPS-PG Ultrapure (LPS from <i>P. gingivalis</i>)	tlr1-pglps
LTA-BS (Standard LTA from <i>B. subtilis</i>)	tlr1-lta

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