# Lipid A from E. coli, Serotype R515 (Re) (TLRGRADE®) (Ready-to-Use)

TLR4 activa View Online » Citations: 8

## **Ordering Information**

**Order Online** »

ALX-581-200-L001	1ml
ALX-581-200-L002	2ml

Manuals, SDS & CofA

**View Online »** 

### **Handling & Storage**

**Use/Stability** As indicated on product label or CoA when stored as recommended.

Handling Keep sterile.

Long Term Storage +4°C

**Shipping** Ambient Temperature

## Regulatory Status RUO - Research Use Only

#### **Product Details**

Activity Activates Toll-like receptor (TLR) 4. Does not activate

TLR2 or other TLRs as determined with splenocytes and

macrophages from TLR4 deficient mice.

Couple Target TLR, TLR4

Couple Type Activator, Ligand

**Formulation** Liquid. Sterile, *ready-to-use* solution in pyrogen-free

double distilled water.

**Purity** ≥99.9% (regarding protein or DNA contaminants; KDO not

detectable)

**Source** Isolated from *E. coli* R515 (Re mutant). Prepared by mild

acid hydrolysis from purified *Escherichia coli* R515 (Re) LPS (isolated by a modification of the PCP extraction method), converted to the uniform salt form and dissolved

in sterile pyrogen-free double distilled water.

Last modified: May 29, 2024

