# LPS from E. coli, Serotype EH100 (Ra) (TLRGRADE®) (Ready-to-Use)

TIPORAPERIAS is commonly used to active TLR4, trigger inflammatory signaling, and co-stimulate the immune response, both *in vitro* and *in vivo*.

Citations: 61

View Online »

# **Ordering Information**

**Order Online** »

ALX-581-010-L001	1ml
ALX-581-010-L002	2ml

Manuals, SDS & CofA

View Online »

- · Potent activator of TLR 4
- TLR4-specific activation as shown in TLR4-deficient cells
- Convenient and stable ready-touse solution
- · Highly cited

## **Handling & Storage**

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

1 year after receipt when stored at +4°C.

Handling Do not ingest. Wear gloves and mask when handling this product! Avoid contact through

all modes of exposure. LPS compounds are highly pyrogenic. Avoid accidental injection; extreme care should be taken when handling in conjunction with hypodermic syringes.

Use must be restricted to qualified personnel.

Long Term Storage +4°C

**Shipping** Ambient Temperature

## Regulatory Status RUO - Research Use Only

### **Product Details**

Alternative Name Lipopolysaccharide from E. coli, Serotype EH100(Ra)

**Appearance** Clear colorless liquid.

Couple Target TLR, TLR4

Couple Type Activator, Ligand

**Formulation** Liquid. Sterile, *ready-to-use* solution in double distilled,

pyrogen-free water.

Purity Absence of detectable protein or DNA contaminants with

agonistic TLR activity.

**Source** Rough (R)-form LPS isolated and purified from *E. coli* 

EH100 (Ra-mutant) by a modification of the PCP

extraction method, converted to the uniform sodium salt form and dissolved in sterile pyrogen-free double distilled

water.

Last modified: May 29, 2024

