## LPS from Salmonella minnesota R345 (Rb) (TLRGRADE<sup>®</sup>) (Ready-to-Use)

TLR4 activa
Citations: 4

View Online »

**Ordering Information** 

Order Online »

ALX-581-015-L002

2ml

Manuals, SDS & CofA

**View Online »** 

## **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**

Activity Strong activator of Toll-like receptor 4 (TLR4). Does not

activate TLR2 or other TLRs as determined with

splenocytes and macrophages from TLR4 deficient mice. No further re-extraction required.Smooth (S)-form LPS are commonly the preferred choice for whole animal studies, whereas Rough (R)-form LPS are primarily used in cellular

in vitro activation studies.

Alternative Name Lipopolysaccharide from Salmonella minnesota R345 (Rb)

Couple Target TLR, TLR4

Couple Type Activator, Ligand

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

double distilled water.

**Purity** Absence of detectable protein or DNA contaminants with

agonistic TLR activity.

**Source** Rough (R)-form LPS, isolated and purified from

Salmonella minnesota R345 (Rb 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



info-

eu@enzolifesciences.com