## LPS from Salmonella minnesota S-Form (TLRGRADE®) (Ready-to-Use)

TLR4 activa
Citations: 6

**Ordering Information** 

Order Online »

ALX-581-020-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 Specific activator of Toll-like receptor 4 (TLR4). Does not

activate TLR2 or other TLRs as determined with

splenocytes and macrophages from TLR4 deficient mice. 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 S-Form

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** Smooth (S)-form LPS isolated and purified from

Salmonella minnesota (wild type) by the hot phenol-water method and dissolved in sterile pyrogen-free double

distilled water.



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989

eu@enzolifesciences.com

Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com