

# LPS from Salmonella abortus equi S-form (TLRGRADE®) (Ready-to-Use) (Biotin)

Useful reagent for evaluating LPS binding and cell activation in white blood cells.

TLR4 activation

Citations: 5

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-581-150-R500

500µl

## Manuals, SDS & CofA

[View Online »](#)

# Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
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. After opening, prepare aliquots and store at -20°C.
Long Term Storage	+4°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Lipopolysaccharide from Salmonella abortus equi S-form
Couple Target	TLR, TLR4
Couple Type	Activator, Ligand
Formulation	Liquid. Sterile, <i>ready-to-use</i> 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 <i>Salmonella abortus equi</i> by modification of the phenol water extraction and PCP method. LPS-Biotin was prepared using the biotin reagent biotinamidocaproate N-hydroxysuccinimide ester. Briefly, LPS at 10mg/ml in distilled water was mixed with biotin reagent in sodium bicarbonate buffer. The reaction mixture was stirred, dialyzed extensively against distilled water in the dark, and sterile filtered.

**Note:** As the biotinylation reagent and standard biotinylation protocol favors protein/peptide (contaminants) about 1000-fold over sugar as a target, it is mandatory to use >99.9% pure LPS to exclude the biotinylation of contaminants. These are usually present in many commercial LPS preparations and would lead to misleading observations of the tracking/staining with the reagents. TLRGRADE<sup>®</sup> LPS fulfills such high quality requirements and the biotinylation procedure is chosen to preserve TLR4 (Toll-like receptor 4) binding and cellular activation using TLRGRADE<sup>®</sup> LPS-Biotin.

For the unconjugated LPS, please see ALX-581-009).

**Last modified: May 29, 2024**



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)