

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

TLR4 activator

Citations: 4

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-581-019-L002	2ml
------------------	-----

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored at +4°C.
<b>Handling</b>	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.
<b>Long Term Storage</b>	+4°C
<b>Shipping</b>	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Activity</b>	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.
<b>Alternative Name</b>	Lipopolysaccharide from <i>Salmonella enteritidis</i> S-Form
<b>Couple Target</b>	TLR, TLR4
<b>Couple Type</b>	Activator, Ligand
<b>Formulation</b>	Liquid. Sterile, <i>ready-to-use</i> solution in pyrogen-free double distilled water.
<b>Purity</b>	Absence of detectable protein or DNA contaminants with agonistic TLR activity.
<b>Source</b>	Smooth (S)-form LPS isolated and purified from <i>Salmonella enteritidis</i> (wild type) by the hot phenol-water method and dissolved in sterile pyrogen-free double distilled water.



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-  
fr@enzolifesciences.com](mailto:info-fr@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)