

# Monophosphoryl Lipid A from Salmonella minnesota R595 (Re) (TLRGRADE<sup>®</sup>) (Ready-to-Use)

Adjuvant

Citations & Value

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-581-202-L001	1ml
ALX-581-2020	500µl

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Keep sterile.
Long Term Storage	+4°C
Shipping	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

Activity	Activates Toll-like receptor (TLR) 4. Does not activate TLR2 as determined with splenocytes and macrophages from TLR4 deficient mice. Exhibits reduced endotoxic and pyrogenic activity compared to lipid A and LPS.
Alternative Name	MPL-A from <i>Salmonella minnesota</i> R595 (Re)
Couple Target	TLR, TLR4
Couple Type	Activator, Ligand
Formulation	Liquid. Sterile, <i>ready-to-use</i> solution in pyrogen-free double distilled water.
Purity	≥99.9% (regarding protein or DNA contaminants; KDO not detectable)
Source	Isolated from <i>Salmonella minnesota</i> R595 (Re mutant). Prepared by acid hydrolysis from purified <i>Salmonella minnesota</i> R595 (Re) LPS (isolated by a modification of the PCP extraction method), converted to the uniform salt form and dissolved in sterile pyrogen-free double distilled water.

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)