

Lipid A from Salmonella minnesota R595 (Re) (TLRGRADE ®) (Ready-to-Use)

TLR4 activa

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-581-201-L002	2ml
------------------	-----

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.
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 mild 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-
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)