

Monophosphoryl Lipid A from E. coli, Serotype R515 (Re) (TLRGRADE®) (Ready-to-Use)

Citations: 3

[View Online »](#)

TLR4 activator

Ordering Information

[Order Online »](#)

ALX-581-203-L001	1ml
ALX-581-203-R500	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 TLR4 (Toll-like receptor 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>E. coli</i> , Serotype R515 (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>E. coli</i> , Serotype R515 (Re mutant). Prepared by acid hydrolysis from purified <i>E. coli</i> , Serotype R515 (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.



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)