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

Citations:16/a

Ordering Information

Order Online »

ALX-581-202-L001	1ml
ALX-581-202-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 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 Salmonella minnesota R595 (Re)

Couple Target TLR, TLR4

Couple Type Activator, Ligand

Formulation Liquid. Sterile, *ready-to-use* solution in pyrogen-free

double distilled water.

Purity ≥99.9% (regarding protein or DNA contaminants; KDO not

detectable)

Source Isolated from *Salmonella minnesota* R595 (Re mutant).

Prepared by acid hydrolysis from purified *Salmonella minnesota* 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 infousa@enzolifesciences.com

European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com