

LPS from E. coli, Serotype O55:B5 (TLRGRADE®) (Ready-to-Use)

TLRGRADE® LPS is commonly used to activate TLR4, trigger inflammatory signaling, and co-stimulate the immune response, both *in vitro* and *in vivo*.

Citations: 74

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-581-013-L001	1ml
ALX-581-013-L002	2ml

Manuals, SDS & CofA

[View Online »](#)

- Potent activator of TLR 4
- TLR4-specific activation as shown in TLR4-deficient cells
- Convenient and stable ready-to-use solution
- Highly cited

Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Lipopolysaccharide from <i>E. coli</i> , Serotype O55:B5 S-form
Couple Target	TLR, TLR4
Couple Type	Activator, Ligand
Formulation	Liquid. Sterile, <i>ready-to-use</i> solution in pyrogen-free double distilled water.
Purity	Absence of detectable protein or DNA contaminants with agonistic TLR activity.
Source	Smooth (S)-form LPS isolated and purified from <i>E. coli</i> O55:B5 by a modification of the phenol water extraction and PCP extraction method, converted to the uniform sodium 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)