

LPS from Salmonella typhimurium S-form (TLRGRADE®) (Ready-to-Use)

TLR4 activat

Citations: 15

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-581-011-L001	1ml
ALX-581-011-L002	2ml

Manuals, SDS & CofA

[View Online »](#)

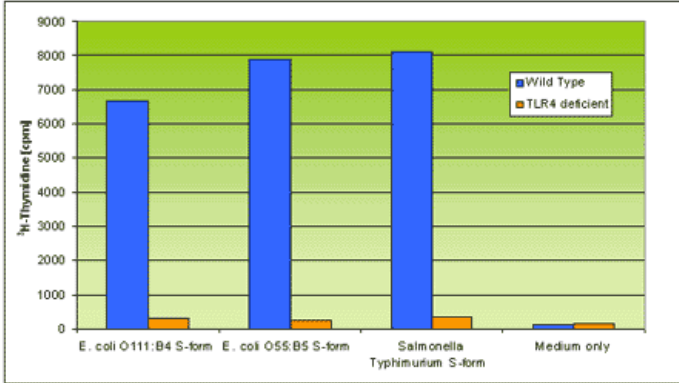


Figure: TLR4-dependent proliferation of mouse splenocytes
Method: 4×10^5 mouse spleen cells were incubated in 0.2ml per well with 10 μ g/ml LPS or with medium only. Proliferation was determined in triplicates by ^3H -thymidine incorporation.

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

Activity	Specific activator of Toll-like receptor (TLR) 4. Does not activate TLR2 or other TLRs as determined with splenocytes and macrophages from TLR4 deficient mice. Smooth (S)-form LPS are commonly the preferred choice for whole animal studies, whereas Rough (R)-form LPS are primarily used in cellular in vitro activation studies.
Alternative Name	Lipopolysaccharide from <i>Salmonella typhimurium</i> S-form
Couple Target	TLR, TLR4
Couple Type	Activator, Ligand
Formulation	Liquid. Sterile, ready-to-use 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>Salmonella typhimurium</i> (wild type) by the hot phenol-water method and dissolved in sterile pyrogen-free double distilled water.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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