

MALP-2 (BULK)

Activator of TLR2/TLR6 complex.

MALP-2 was originally isolated from *Mycoplasma fermentans*. This MALP-2 corresponds to the originally isolated isomer, which expresses potent endotoxin-like activity and approaches in certain experimental systems the toxicity of LPS. For description of the stereochemistry of MALP-2. MALP-2 signalling, unlike that of LPS, is not transduced via TLR4, but is induced via TLR2 and TLR6 signalling. For more information about MALP-2 see www.malp-research.de. BULK size (500µg) of biologically active, pyrogen-free and sterile MALP-2 for *in vitro* and animal *in vivo* studies.

Citations: 13

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-162-034-C500	500µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 3 years when stored at -20°C. This stock solution is stable for many months when kept in the refrigerator. MALP-2 solutions in the presence of carrier (50mM octyl glucoside, see Background/Technical Information below) are stable at +4°C at least 6 weeks. Some (10%) decrease of activity may occur shortly after transfer into new vessels. This is not due to decomposition but to adsorption (see below). Storage at room temperature for short periods of time does not influence bioactivity. When deep-frozen, MALP-2 does not lose activity within years.
Handling	Avoid accidental injection, extreme care should be taken with hypodermic syringes. Avoid inhalation and prevent this compound from entering the bloodstream. Avoid freeze/thaw cycles.
Short Term Storage	-20°C
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Macrophage-activating lipopeptide-2, S-[2,3-bis(Palmitoyloxy)-(2R)-propyl-cysteiny]-GNNDESNISFKEK]
Concentration	1mg/ml
Formula	$C_{99}H_{167}N_{19}O_{30}S$
Formulation	Liquid. Pyrogen-free, sterile solution in 33% (v/v) 2-propanol/water.
MW	2135.2
Purity Detail	HPLC purified
Source	Synthetic.

MALP-2 sticks avidly to glass or plastic. Since it is active at very low concentrations, i.e. at high dilutions, these low amounts of material tend to become adsorbed to pipette tips and plastic or glass containers. To avoid this, prepare several dilution steps of the MALP-2 stock solution with medium containing 5% autologous serum or buffers with 2% HSA. The presence of 50mM octyl glucoside in the first dilution step is often beneficial and gives rise to higher activity. Avoid small volumes in large vessels. Do not filter. Because of the ester-bound fatty acids, MALP-2 is sensitive to alkali or acid. The medium should be preconditioned in the CO₂ incubator in order to ensure that the pH does not become alkaline. Add cells in preconditioned medium as soon as possible and incubate.

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-
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)