

Ac-muramyl-Ala-D-Glu-amide

NALP3 ligand

Specific ligand for NALP3. For inactive control compound see MDP-DD (Prod. No. ALX-151-036).

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	MDP-LD, N-Acetylmuramyl-L-alanyl-D-isoglutamine, Adjuvant peptide
Appearance	White to off-white powder.
CAS	53678-77-6
Formula	$C_{19}H_{32}N_4O_{11}$
MW	492.5
Purity	≥95% (HPLC)
Solubility	Soluble in water (50mg/ml).
Source	Synthetic.

Last modified: May 29, 2024



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