

MDP

Muramyl dipeptide, L-D isomer; NOD2 ligand

Catalog code: tlrl-mdp

<https://www.invivogen.com/mdp>

For research use only

Version 23G26-MM

PRODUCT INFORMATION

Contents

- 5 mg Muramyl dipeptide (MDP)
- 1.5 ml endotoxin-free water

Storage and stability

- MDP is shipped at room temperature. Upon receipt, store at -20°C.
- Upon resuspension, prepare aliquots of MDP and store at -20°C. Resuspended product is stable for 6 months at -20°C when properly stored. Avoid repeated freeze-thaw cycles.

Quality control

- Activation of NOD2 has been confirmed using HEK-Blue™ NOD2 cells.
- The absence of NOD1 activity has been confirmed using HEK-Blue™ NOD1 cells.
- The absence of bacterial contamination (e.g. lipoproteins and endotoxins) has been confirmed using HEK-Blue™ TLR2 and HEK-Blue™ TLR4 cells.

DESCRIPTION

Muramyl dipeptide (MDP) is a synthetic immunoreactive peptide composed of N-acetylmuramic acid linked by its lactic acid moiety to the N-terminus of an L-alanine D-isoglutamine dipeptide. MDP is the minimal bioactive peptidoglycan motif found in almost all bacteria. It was first identified as an active component in Freund's complete adjuvant¹.

MDP is recognized by the cytosolic receptor NOD2^{2,3}. Ligand-bound NOD2 oligomerizes and signals via the serine/threonine RIP2 (RICK,CARDIAK) kinase through CARD-CARD homophilic interactions⁴. Once activated, RIP2 mediates ubiquitination of NEMO/IKK γ leading to the activation of NF- κ B and the production of inflammatory cytokines. Furthermore, poly-ubiquitinated RIP2 recruits TAK1, which leads to IKK complex activation and the activation of MAPKs⁵. This signaling involves the adaptor protein CARD9⁶.

NOD2 recognition of MDP is stereospecific to the L-D isomer, excluding any reaction to D-D or L-L analogs³. The potent adjuvant activity of MDP may also be linked to an activation of the NLRP3 inflammasome⁷. Of note, NOD2 mutants associated with susceptibility to Crohn's disease have been found to be deficient in their recognition of MDP^{2,3}.

1. Ogawa C. *et al.*, 2011. Muramyl dipeptide and its derivatives: peptide adjuvant in immunological disorders and cancer therapy. *Curr Bioact Compd.* 7(3):180-97. 2. Girardin S.E. *et al.*, 2003. Nod2 is a general sensor of peptidoglycan through muramyl dipeptide (MDP) detection. *J Biol Chem.* 278(11):8869-72. 3. Inohara N. *et al.*, 2003. Host recognition of bacterial muramyl dipeptide mediated through NOD2. Implications for Crohn's disease. *J Biol Chem.* 278(8):5509-12. 4. Kobayash K. *et al.*, 2002. RICK/Rip2/CARDIAK mediates signalling for receptors of the innate and adaptive immune systems. *Nature* 416:194-9. 5. Kobayashi K. *et al.*, 2005. Nod2-dependent regulation of innate and adaptive immunity in the intestinal tract. *Science* 307: 731-734. 6. Hsu Y. *et al.*, 2007. The adaptor protein CARD9 is required for innate immune responses to intracellular pathogens. *Nat Immunol.* 8(2):198-205. 7. Martinon F. *et al.*, 2004. Identification of bacterial muramyl dipeptide as activator of the NALP3/cryopyrin inflammasome. *Curr Biol.* 14(21):1929-34.

TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Asia: +852 3622-3480

E-mail: info@invivogen.com

PRODUCT PROPERTIES

Synonym: N-Acetylmuramyl-L-Alanyl-D-Isoglutamine

CAS number: 53678-77-6

Formula: C₁₉H₃₂N₄O₁₁

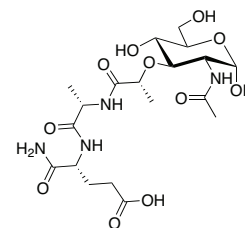
Molecular weight: 492.5 g/mol

Specificity: NOD2

Solubility: 10 mg/ml in water

Working concentration: 10 ng-10 μ g/ml

Chemical structure:



METHODS

Preparation of stock solution (10 mg/ml)

1. Add 500 μ l endotoxin-free water (provided) to the vial containing 5 mg of MDP and vortex to solubilize.

NOD2 activation using MDP

MDP can be used to activate NOD2 in cells expressing this receptor, such as HEK-Blue™ NOD2 cells. These cells express the human or murine NOD2 gene and an NF- κ B inducible SEAP reporter gene. Levels of SEAP can be easily determined using HEK-Blue™ Detection, a cell culture medium that allows the detection of SEAP as the reporter protein is secreted by the cells. For more information, visit: <https://www.invivogen.com/hek-blue-nod>.

1. Dispense 20 μ l of MDP at various concentrations (10 ng-10 μ g/ml) per well of a 96-well plate.
2. Prepare a cell suspension ~280,000 cells per ml in HEK-Blue™ Detection.
3. Add 180 μ l of the cell suspension (~50,000 cells) to each MDP-containing well.
4. Incubate the plate for 6-24 h at 37°C, 5% CO₂.
5. Determine SEAP levels using a spectrophotometer at 620-655 nm.

RELATED PRODUCTS

Product	Description	Cat. Code
HEK-Blue™ Detection	SEAP Detection reagent	hb-det2
HEK-Blue™ hNOD2 Cells	Human NOD2 reporter cells	hkb-hnod2
HEK-Blue™ mNOD2 Cells	Murine NOD2 reporter cells	hkb-mnod2
MDP Control	L-L isomer, negative control	tlrl-mdpctl