

# C12-iE-DAP

NOD1 ligand

Catalog code: tlr1-c12dap

<https://www.invivogen.com/c12-ie-dap>

For research use only

Version 24C11-MM

## PRODUCT INFORMATION

### Contents

- 1 mg C12-iE-DAP.

*Note:* It is a mixture of Lauroyl- $\gamma$ -D-Glu-D-mDAP and Lauroyl- $\gamma$ -D-Glu-L-mDAP.

- 1.5 ml endotoxin-free water

### Storage and stability

- C12-iE-DAP is provided lyophilized and shipped at room temperature. Upon receipt, store at -20°C.

- Upon resuspension, prepare aliquots of C12-iE-DAP and store at -20°C. Resuspended product is stable for 1 year when properly stored. Avoid repeated freeze-thaw cycles.

### Quality control

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

## DESCRIPTION

C12-iE-DAP is a chemically synthesized NOD1 agonist. It is an acylated derivative of the dipeptide iE-DAP ( $\gamma$ -D-Glu-mDAP), present in the peptidoglycan of a subset of bacteria that include Gram-negative bacilli and some Gram-positive bacteria such as *Bacillus subtilis* and *Listeria monocytogenes*<sup>1</sup>. C12-iE-DAP was generated by the addition of a lauroyl (C12) group to the glutamic residue of iE-DAP. Similar to iE-DAP, C12-iE-DAP is recognized by NOD1 (CARD4), an intracellular sensor expressed in multiple tissues, including intestinal epithelia cells. Recognition of C12-iE-DAP by NOD1 induces a signaling cascade involving the serine/threonine RIP2 (RICK, CARDIAK) kinase, which interacts with IKK to trigger the activation of NF- $\kappa$ B and the production of inflammatory cytokines such as TNF- $\alpha$  and IL-6<sup>2</sup>. C12-iE-DAP stimulates NOD1 at concentrations 100- to 1000-fold lower than does iE-DAP.

1. Chamailard M. *et al.*, 2003. An essential role for NOD1 in host recognition of bacterial peptidoglycan containing diaminopimelic acid. *Nat. Immunol.*4(7):702-7.
2. Park JH. *et al.*, 2007. RICK/RIP2 mediates innate immune responses induced through Nod1 and Nod2 but not TLRs. *J Immunol.* 178(4):2380-6.

## CHEMICAL PROPERTIES

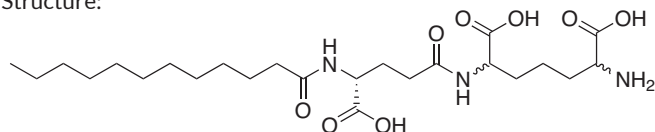
**Synonym:** Lauroyl- $\gamma$ -D-glutamyl-meso-diaminopimelic acid

**Formula:** C<sub>24</sub>H<sub>43</sub>N<sub>3</sub>O<sub>8</sub>

**Molecular weight:** 501.61 g/mol

**Solubility:** 10 mg/ml DMSO or methanol

**Structure:**



## METHOD

### Preparation of stock solution (1 mg/ml)

1. Add 100  $\mu$ l DMSO (or methanol) to 1 mg of C12-iE-DAP.
2. Vortex until completely resuspended.
3. Once C12-iE-DAP is resuspended, add 900  $\mu$ l endotoxin-free water (provided).
4. Prepare aliquots and store at -20°C. Further dilutions can be prepared by adding the appropriate amount of endotoxin-free water.

### NOD1 activation using C12-iE-DAP

C12-iE-DAP can be used to activate NOD1 in cells expressing this receptor such as HEK-Blue™ NOD1 cells. These cells were designed to study NOD1 stimulation by monitoring NF- $\kappa$ B activation. Stimulation of HEK-Blue™ NOD1 cells with a NOD1 agonist activates NF- $\kappa$ B, which induces the production of SEAP (secreted embryonic alkaline phosphatase). Levels of SEAP can be easily determined using HEK-Blue™ Detection, a cell culture medium that allows the detection of SEAP as it is secreted by the cells. For more information visit: <https://www.invivogen.com/hek-blue-nod1>.

1. Add 20  $\mu$ l of C12-iE-DAP at various concentrations (10 ng-10  $\mu$ g/ml) per well of a 96-well plate.
2. Prepare a cell suspension (~280,000 cells per ml) in HEK-Blue™ Detection medium and immediately add 180  $\mu$ l of the cell suspension (~50,000 cells) to each C12-iE-DAP-containing well.
3. Incubate the plate for 6-24 h at 37°C, 5% CO<sub>2</sub>.
4. Determine SEAP levels using a spectrophotometer at 620-655 nm.

## RELATED PRODUCTS

Product	Description	Cat.Code
HEK-Blue™ hNOD1 cells	Human NOD1 reporter cells	hkb-hnod1
HEK-Blue™ mNOD1 cells	Murine NOD1 reporter cells	hkb-mnod1
HEK-Blue™ Detection	SEAP detection medium	hb-det2
iE-DAP	NOD1 ligand	tlr1-dap
Tri-DAP	NOD1 ligand	tlr1-tdap

## 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](mailto:info@invivogen.com)