



## Product Data Sheet

---

**Product Name:** Ala- $\gamma$ -D-Glu-DAP

**Catalog Number:** AS-60774 (1 mg)                      **Lot Number:** See label on vial

**Sequence:** H-Ala-D- $\gamma$ -Glu-diaminopimelic acid (3-letter code)  
A-( $\gamma$ -e)-DAP (1-letter code)

**Molecular Weight:** 391.4

**% Peak Area by HPLC:**  $\geq 95$

**Appearance:** Lyophilized white solid

**Peptide Reconstitution:** Ala- $\gamma$ -D-Glu-DAP peptide is freely soluble in H<sub>2</sub>O.

**Storage:** Ala- $\gamma$ -D-Glu-DAP peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at  $-20^{\circ}\text{C}$  or lower. Reconstituted peptide can be aliquoted and stored at  $-20^{\circ}\text{C}$  or lower.

**Description:** This peptide, with the diaminopimelic acid coupled to the  $\gamma$ -carboxylic acid of the D-isomer of Glu, stimulates Nod1-dependent apoptosis. Ref: da Silva Correia, J. et al. *Proc. Natl. Acad. Sci. USA* **103**, 1840 (2006); Baum, EZ. et al. *Antimicrobial Agents Chemo.* **50**, 230 (2006).

**Additional Information:** Listed below are relevant information that may provide a guideline on how to use this product. End users will have to adapt to their own specific applications.

$\gamma$ TriDAP (Ala- $\gamma$ Glu-mesoDAP) was chemically synthesized by Anaspec (San Jose, CA). Cells were left untreated or treated with  $\gamma$ TriDAP or MDP (20  $\mu\text{g}/\text{ml}$  each) in the presence or absence of IFN- $\gamma$  (1,000 units/ml) for 24 h-[Da Silva Correia, J. et al. Proc. Natl. Acad. Sci. 104, 6764 \(2007\).](#)

**Published Citations:**

Da Silva Correia, J. et al. *Proc. Natl. Acad. Sci.* **104**, 6764 (2007).

*For Research Use Only*