

## **Product Data Sheet**

Product Name: GALA, Pore-Forming Peptide

Catalog Number: AS-62311 (1 mg) Lot Number: See label on vial

Sequence: H-Trp-Glu-Ala-Ala-Leu-Ala-Glu-Ala-Leu-Ala-Glu-Ala-Leu-Ala-Glu-His-

Leu-Ala-Glu-Ala-Leu-Ala-Glu-Ala-Leu-Glu-Ala-Leu-Ala-Ala-OH

(3-letter code)

WEAALAEALAEALAEHLAEALAEALEALAA (1-letter code)

Molecular Weight: 3032.4

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: GALA, Pore-Forming peptide is freely soluble in 1% NH<sub>4</sub>OH.

Storage: GALA, Pore-Forming peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at –20°C or lower. Reconstituted peptide can be aliquoted and stored at – 20°C or lower.

Description: GALA is a 30 amino acid synthetic peptide with a glutamic acid-alanine-leucine-alanine (EALA) repeat. It also contains a histidine and tryptophan residue as spectroscopic probes. This peptide was designed to explore how viral fusion protein sequences interact with membranes. It was used to study the importance of the helix length, hydrophobicity, and hydrophobic moment on the formation, structure, and function of ion channels. The membrane-interacting properties of GALA have been extensively documented. Investigations with analogs of cytotoxic peptides clarify the importance of peptide charge and the role of particular amino acids or arrays of amino acids in the conformation, membrane-binding affinity, and pore-forming abilities of these toxins. Ref: Francois, N. et al. Biophys. J. 76, 2121 (1999), Li, W. et al. *Adv. Drug Del. Rev.* 56, 967 (2004). Ref: Francois, N. et al. *Biophys. J.* 76, 2121 (1999), Li, W. et al. *Adv. Drug Del. Rev.* 56, 967 (2004).

## Related Products:

NameCat #SizeApidaecin IBAS-62044 1 mg

**GNNRPVYIPQPRPPHPRL** 

Magainin 2 AS-20639 0.5 mg
GIGKFLHSAKKFGKAFVGEIMNS AS-20640 1 mg