

Product Data Sheet

Product Name: Magainin 2

Catalog Number: AS-20639 (0.5 mg) Lot Number: See label on vial

AS-20640 (1 mg)

Sequence: H-Gly-Ile-Gly-Lys-Phe-Leu-His-Ser-Ala-Lys-Lys-Phe-Gly-Lys-Ala-Phe-

Val-Gly-Glu-Ile-Met-Asn-Ser-OH (3-letter code) GIGKFLHSAKKFGKAFVGEIMNS (1-letter code)

Molecular Weight: 2467.9

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Using H2O, reconstitute by adding 50-70 µl to 0.5 mg Magainin 2

peptide.

Storage: Magainin 2 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: Magainin 2 assumes an amphiphilic helix when bound to acidic phospholipids, forming a pore composed of a dynamic, peptide-lipid supramolecular complex. Ref: Zasloff, M. *Proc. Natl. Acad. Sci. USA.* **84**, 5449 (1987); Cruciani, RA. et al. *EJPMOL* **8**, 187 (1992); Corzo, G. et al. *Biochem. J.* **359**, 35 (2001); Matsuzaki, K. et al. *Biochem.* **36**, 2104 (1997).

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.

The LL-37 and magainin 2 peptides were obtained from Anaspec. Using the *Escherichia coli* K12 JM109 strain (New England Biolabs), we measured MIC values of 3 and 12 μ M for LL-37 and magainin 2, respectively. These values were consistent with those measured previously for these peptides with other *E. coli* strains . For the cell viability/permeabilization assays, cultures of *E. coli* JM109 were grown exponentially in Luria–Bertani (LB) medium (OD_{600nm} = 0.6) and then pelleted by centrifugation at 1600g for 10 min at 4 °C-Bourbon, C. et al. *Anal. Biochem.* **381**, 279 (2008).

Published Citations:

Maher, S. et al. *Biochem. Pharmacol.* **71**, 1289 (2006). Bourbon, C. et al. *Anal. Biochem.* **381**, 279 (2008). Palmero, EF, et al. *Biomacromolecules* **10**, 1416 (2009).

Related Products:

 Name
 Cat #
 Size

 Magainin 1
 AS-20791
 0.5 mg

 (GIGKFLHSAGKFGKAFVGEIMKS)
 AS-20792
 1 mg

For Research Use Only