

Product Name:	Magainin 2	
Catalog Number:	AS-20639 (0.5 mg) AS-20640 (1 mg)	Lot Number: See label on vial
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 5	50-70 μl to 0.5 mg Magainin 2

peptide Reconstitution. Using H2O, reconstitute by adding 50-70 µr 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  $\mu$ 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<sub>600nm</sub> = 0.6) and then pelleted by centrifugation at 1600*g* 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 # S	ize
Magainin 1 (GIGKFLHSAGKFGKAFVGEIMKS)	AS-20791 AS-20792	0.5 mg 1 mg <i>For Research Use Only</i>