



## Product Data Sheet

---

**Product Name:** Melittin, honey bee

**Catalog Number:** AS-62366 (1 mg)                      **Lot Number:** See label on vial

**Sequence:** H-Gly-Ile-Gly-Ala-Val-Leu-Lys-Val-Leu-Thr-Thr-Gly-Leu-Pro-Ala-Leu-Ile-Ser-Trp-Ile-Lys-Arg-Lys-Arg-Gln-Gln-OH (3-letter code)  
GIGAVLKVLTTGLPALISWIKRKRQQ (1-letter code)

**Molecular Weight:** 2848.5

**% Peak Area by HPLC:** ≥ 95

**Appearance:** Lyophilized white powder

**Peptide Reconstitution:** Using H<sub>2</sub>O, reconstitute by adding > 100 µl to 1 mg Melittin peptide.

**Storage:** Melittin 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:** Melittin, a 26-residue bee venom peptide, is known to induce murine antibodies specific for its hydrophilic C-terminus of residues 20 to 26 and T-cell responses specific for its hydrophobic mid-region (residue 11 to 19). This peptide is an anti-inflammatory agent, it inhibits the Lyme disease spirochete. Smith, R. et al. *J. Mol. Biol.* **241**, 456 (1994), King, T. et al. *J. Immunol.* **153**, 1124 (1994).

### Related Products:

Name	Cat #	Size
Apamin (CNCKAPETALCARRCQQH-NH <sub>2</sub> (Disulfide bridge: 1-11; 3-15))	AS-60772	0.5 mg

*For Research Use Only*