

Product Data Sheet

Product Name: Cecropin A

Catalog Number: AS-24009 (0.5 mg) Lot Number: See label on vial

AS-24010 (1 mg)

Sequence: H-Lys-Trp-Lys-Leu-Phe-Lys-Lys-Ile-Glu-Lys-Val-Gly-Gln-Asn-Ile-Arg-

Asp-Gly-Ile-Ile-Lys-Ala-Gly-Pro-Ala-Val-Ala-Val-Val-Gly-Gln-Ala-Thr-Gln-

Ile-Ala-Lys-NH2 (3-letter code)

KWKLFKKIEKVGQNIRDGIIKAGPAVAVVGQATQIAK-NH2 (1-letter

code)

Molecular Weight: 4004.8 % Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Cecropin A peptide is freely soluble in H₂O.

Storage: Cecropin A 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: Cecropin A is a naturally occurring, linear, cationic, 37-residue antimicrobial peptide. Cecropin A kills bacteria by dissipating transmembrane electrochemical ion-gradients.

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.

Antimicrobial peptides cecropin A and B and magainin-1 were received from Anaspec (San Jose, CA). Capture solutions containing AMPs or antibodies diluted in PBS were injected into the PDMS channels and incubated overnight at 4°C. Concentrations of the recognition molecules in the capture solutions were as follows: Polymyxins B and E, 10 mg/ml; cecropins A, B and P, magainin-1, and parasin, 1 mg/ml; melittin and bactenecin, 250 µg/ml; control and target-specific antibodies, 10 µg/ml. After overnight incubation with the capture solutions, the channels were emptied and rinsed with PBS. The slides were then blocked for 30 min in 10 mg/ml gelatin in PBS, dried, and stored at 4°C for up to 2 weeks-Kulagina, NV. et al. Sens Actuators B Chem. 121, 150 (2007).

Published Citations:

Kulagina, NV. et al. *Analytica. Chimica. Acta.* **575**, 9 (2006) Kulagina, NV. et al. *Sens Actuators B Chem.* **121**, 150 (2007).

Related Products:

 Name
 Cat # Size

 Cecropin B
 AS-24011 0.5 mg

(KWKVFKKIEKMGRNIRNGIVKAGPAIAVLGEAKAL-NH2)

For Research Use Only