

## **Product Data Sheet**

Product Name: mCRAMP, mouse

Catalog Number: AS-61305 (1 mg) Lot Number: See label on vial

Sequence: H-Gly-Leu-Leu-Arg-Lys-Gly-Glu-Lys-Ile-Gly-Glu-Lys-Leu-Lys-Lys-Ile-

Gly-Gln-Lys-Ile-Lys-Asn-Phe-Phe-Gln-Lys-Leu-Val-Pro-Gln-Pro-Glu-Gln-

OH (3-letter code)

GLLRKGGEKIGEKLKKIGQKIKNFFQKLVPQPEQ (1-letter code)

Molecular Weight: 3879.7

% 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 mCRAMP peptide.

Storage: mCRAMP 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: This cathelicidin-related antimicrobial peptide (mCRAMP) is the sole murine cathelicidin. mCRAMP expression in the intestinal tract is restricted to surface epithelial cells in the colon. mCRAMP shows antimicrobial activity against the murine enteric pathogen *Citrobacter rodentium* and destroys skin *Candida albicans*. Limura, M. et al. *J. Immunol.* **174**, 4901 (2005); Lopez-Garcia, B. et al. *J. Invest. Dermatol.* **125**, 108 (2005).

## **Published Citations:**

Gryllos, I. et al. PNAS 105, 16755 (2008).

For Research Use Only