



Product Data Sheet

Product Name: MAPS, Core, 8-Branch

Catalog Number: AS-21006 (10 mg) **Lot Number:** See label on vial

Sequence: H-(Lys)₄-(Lys)₂-Lys-Ala-NH₂ (3-letter code)
K₄K₂KA-NH₂ (1-letter code)

Molecular Weight: 985.3

Peptide Purity: >95%

Appearance: Lyophilized white crystal.

Peptide Reconstitution: MAPS peptide is freely soluble in water.

Storage: MAPS 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.

Published Citations:

[Abboud, N. et al. *J Biol Chem* **284**, 25077 \(2009\).](#)

For Research Use Only