

Product Data Sheet

Product Name: Proadrenomedullin N-terminal 20 peptide, PAMP (1-20)

Catalog Number: AS-65455-5 (5 mg) Lot Number: See label on vial

Sequence: H-Ala-Arg-Leu-Asp-Val-Ala-Ser-Glu-Phe-Arg-Lys-Lys-Trp-Asn-Lys-Trp-

Ala-Leu-Ser-Arg-NH2

ARLDVASEFRKKWNKWALSR-NH2

Molecular Weight: 2460.9

Peptide Purity: Peak Area by HPLC ≥95%

Storage: This 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: PAMP (1-20) is a biologically active molecule produced by

posttranslational processing of proadrenomedullin. This peptide is found in a variety of tissues and exerts important biological activities through specific cell surface receptors. PAMP (1-20) is involved in many physiological processes, including vasodilation, bronchodilation, hormone secretion, cell migration, apoptosis, and angiogenesis. This peptide has been implicated in the pathophysiology of several diseases such as renal and cardiovascular disorders, stroke, cancer, and

diabetes.

Related Products:

Adrenomedullin (1 - 52), human, Cat# AS-60447 Adrenomedullin (22 - 52), human, Cat# AS-60449

References:

1. Mahata, M. et al. *Hypertension* **32**, 907 (1998)

2. Kamohara, M. et al. *Biochem Biophys Res Commun* **330**,1146

(2005)

3. López, J. et al. *Int Rev Cytol* **221**,1 (2002)

For Research Use Only