



Product Data Sheet

Product Name: Adrenomedullin (1-50), rat

Catalog Number: AS-60454 (0.5 mg) **Lot Number:** See label on vial

Sequence: H-Tyr-Arg-Gln-Ser-Met-Asn-Gln-Gly-Ser-Arg-Ser-Thr-Gly-Cys-Arg-Phe-Gly-Thr-Cys-Thr-Met-Gln-Lys-Leu-Ala-His-Gln-Ile-Tyr-Gln-Phe-Thr-Asp-Lys-Asp-Lys-Asp-Gly-Met-Ala-Pro-Arg-Asn-Lys-Ile-Ser-Pro-Gln-Gly-Tyr-NH₂ (Disulfide bridge: 14-19) (3-letter code)
YRQSMNQGSRSTGCRFGTCTMQKLAHQIYQFTDKDKDGMAPRNKISP
QGY-NH₂ (Disulfide bridge: 14-19) (1-letter code)

Molecular Weight: 5730.5

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Using H₂O, reconstitute by adding 50 µl to 0.5 mg Adrenomedullin peptide.

Storage: Adrenomedullin 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: Rat adrenomedullin, rADM, (1-50) and its C-terminal rADM (11-50) induce a dose-dependent and endothelium-independent vasodilation on the arterial mesenteric vasculature. Ref: Belloni, AS. et al. *Life Sci.* **63**, 2313 (1998); Sone, M. et al. *Peptides* **18**, 1125 (1997).

Related Products:

Name	Cat #	Size
Adrenomedullin (1-52), human (YRQSMNMFQGLRSFGCRFGTCTVQKLAHQIYQFTDKDKDNVAPRSKISPQGY-NH ₂) (Disulfide bridge: 16-21)	AS-60447	0.5 mg

For Research Use Only