

## **Product Data Sheet**

Product Name: Angiotensin I, human

Catalog Number: AS-20627 (5 mg) Lot Number: See label on vial

AS-20628 (25 mg)

Sequence: H-Asp-Arg-Val-Tyr-lle-His-Pro-Phe-His-Leu-OH (3-letter code)

DRVYIHPFHL (1-letter code)

Molecular Weight: 1296.5 % Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Use distilled or higher quality water. Add water directly to the lyophilized peptide powder to obtain a final concentration of approximately 0.5 mg/mL to 1mg/mL or less. Gently vortex to mix. For peptides that have poor solubility in the suggested solvent, brief sonication may increase solubility in some cases.

Storage: Peptide is shipped at ambient temperature. Upon receipt, store lyophilized powder at – 20°C or lower. Reconstituted peptide should be aliquoted into several freezer vials and stored at –20°C or lower. Do not freeze thaw.

Description: This human Angiotensin I (Ang I) sequence also corresponds to horse, sheep, pig, and rat Ang I. Ang I is cleaved to Ang II by the angiotensin-converting enzyme (ACE). There is also evidence for non-angiotensin-converting enzyme-dependent conversion of Ang I to Ang II. Human chymase efficiently converts the 10-mer Ang I to the 8-mer hormone Ang II by splitting the Phe8-His9 bond in Ang I. Ref: Lundequist, A. et al. *J Biol Chem* **279**, 32339 (2004); Olson, S. et al. *Am J Physiol Lung Cell Mol Physiol* **287**, L559 (2004); Sanker, S. et al. *J Biol Chem* **272**, 2963 (1997).

## **Published Citations:**

Zhang, C. et al. *Clin Exp Pharmacol Physiol* **38**, 55 (2011). Keppel, T. et al. *Rapid Comm Mass Spectrom* **24**, 6 (2009). Rauniyar, N. et al. *Anal Chem* **81**, 782 (2009).

For Research Use Only