



Product Data Sheet

Product Name: Angiotensin II, human
Catalog Number: AS-20633 (5 mg) Lot Number: See label on vial
AS-20634 (25 mg)

Sequence: H-Asp-Arg-Val-Tyr-Ile-His-Pro-Phe-OH (3-letter code)
DRVYIHPF (1-letter code)

Molecular Weight: 1046.2

% 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: The octapeptide angiotensin II (Ang II) exerts a wide range of effects on the cardiovascular system. It is also implicated in the regulation of cell proliferation, fibrosis and apoptosis. Ang II is formed by cleavage of Ang I by the angiotensin-converting enzyme (ACE) or chymases. Human heart chymase, a chymotrypsin-like serine proteinase, hydrolyzes the Phe8-His9 bond to yield the octapeptide hormone angiotensin II and His-Leu. Ref: Lundquist, A. et al. *J Biol Chem* **279**, 32339 (2004); Spyroulias, G. et al. *Eur J Biochem* **270**, 2163 (2003); Cobb, C. et al. *Gen Comp Endocrinol* **127**, 8 (2002); Kinoshita, A. et al. *J Biol Chem* **266**, 19192 (1991)

Published Citations:

- Zhang, C. et al. *Clin Exp Pharmacol Physiol* **38**, 55 (2011).
- Badu-Tawiah, A. et al. *J Am Soc Mass Spectrom* **21**, 1423 (2010).
- Zhao, Y. et al. *J Cell Physiol* 10.1002/jcp.22209 (2010).
- Berkout, V. et al. *Anal Chem* **81**, 725 (2009).
- Kim, J. et al. *J Cerebral Blood Flow Metab* **29**, 621 (2009).
- Rauniyar, N. et al. *Anal Chem* **81**, 782 (2009).
- Gardner, M. et al. *Anal Chem* **80**, 4807 (2008).
- Zhao, Y. et al. *Acta Biochim Biophys Sin* **40**, 304 (2008).
- Chipitsyna, G. *Endocrinology* **148**, 2198 (2007).
- Zhang, S. et al. *J Proteome Res* **6**, 1700 (2007).
- Zhang, S. et al. *J Anal App Pyrolysis* **80**, 353 (2007).
- Crowe, M.C. et al. *JASMS* **15**, 1581 (2004).
- Woods, A. S. et al. *J Biomol Tech* **14**, 1 (2003).

For Research Use Only