



Product Data Sheet

Product Name: BNP-32, human

Catalog Number: AS-24015 (0.5 mg) Lot Number: See label on vial
AS-24016 (1 mg)

Sequence: H-Ser-Pro-Lys-Met-Val-Gln-Gly-Ser-Gly-Cys-Phe-Gly-Arg-Lys-Met-Asp-Arg-Ile-Ser-Ser-Ser-Ser-Gly-Leu-Gly-Cys-Lys-Val-Leu-Arg-Arg-His-OH
(Disulfide bridge: 10-26) (3-letter code)
SPKMQGSGCFGRKMDRISSSSSGLGCKVLRH
(Disulfide bridge: 10-26) (1-letter code)

Molecular Weight: 3465.1

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: BNP-32 is freely soluble in H₂O.

Storage: BNP-32 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: This 32 amino acid peptide contains a 17 amino acid ring structure that is common to all natriuretic peptides. It is also called the brain natriuretic peptide (BNP) because it was first identified in porcine brain; however, the main source of this peptide is not the brain but the cardiac ventricle. This cardiac neurohormone is secreted from the ventricles in response to volume expansion and pressure overload. It has natriuretic and vasodilatory effects and suppresses the renin-angiotensin-aldosterone system. Ref: Cardarelli, R. et al. *J. Am. Board Fam. Prac.* **16**, 327 (2003); Sudoh, T. et al. *Nature* **332**, 78 (1988); Cheung, B. et al. *JAMA* **280**, 1983 (1998); Struthers, A. *BMJ* **308**, 1615 (1994); De Lemos, JA. et al. *Lancet* **362**, 316 (2005); Kambayashi, Y. et al. *Biochem. Biophys. Res. Commun.* **159**, 1427 (1989).

For Research Use Only