



## Product Data Sheet

---

**Product Name:** NGR Peptide 1

**Catalog Number:** AS-60189-1 (1 mg)                      **Lot Number:** See label on vial

**Sequence:** H-Cys-Asn-Gly-Arg-Cys-Gly-Gly-D-Lys-D-Leu-D-Ala-D-Lys-D-Leu-D-Ala-D-Lys-D-Lys-D-Leu-D-Ala-D-Lys-D-Leu-D-Ala-D-Lys-NH<sub>2</sub> (Disulfide bridge: 1-5) (3-letter code)  
CNGRCGGklaklakkklakl-NH<sub>2</sub> (Disulfide bridge: 1-5) (1-letter code)

**Molecular Weight:** 2168.7

**Peptide Purity:** >95%

**Appearance:** Lyophilized white powder

**Peptide Reconstitution:** NGR Peptide is freely soluble in water.

**Storage:** NGR 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.

**References:** Ellerby, HM. et al. *Nature Medicine* **5**, 1032 (1999); Colombo, G. et al. *J. Biol. Chem.* **277**, 47891 (2002); Van Hensbergen, Y. et al. *Biochem. Pharmacol.* **63**, 897 (2002); Plesniak, L. et al. *Protein Sci.* **13**, 1988 (2004).

*For Research Use Only*