



## Product Data Sheet

---

**Product Name:** Atrial Natriuretic Peptide (1 - 28), rat

**Catalog Number:** AS-20651 (0.5 mg)                      Lot Number: See label on vial  
AS-20652 (1 mg)

**Sequence:** H-Ser-Leu-Arg-Arg-Ser-Ser-Cys-Phe-Gly-Gly-Arg-Ile-Asp-Arg-Ile-Gly-  
Ala-Gln-Ser-Gly-Leu-Gly-Cys-Asn-Ser-Phe-Arg-Tyr-OH  
(Disulfide bridge: 7-23) (3-letter code)  
SLRRSSCFGGRIDRIGAQSGLGCNSFRY (Disulfide bridge: 7-23)  
(1-letter code) (1-letter code)

**Molecular Weight:** 3062.5

**% 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:** Atrial natriuretic peptide (ANP 1-28, rat) helps regulate blood pressure and sodium/water balance under normal and pathophysiological environment. Ref: De Paolis, P. et al. Annals Clin Lab Sci 37, 135 (2007).

:

***For Research Use Only***