

Product Name:	C-Type Natriuretic Peptide (32-53), human, porcine		
Catalog Number:	AS-24244 (1 mg)	Lot Number: See label on vial	
Sequence:	H-Gly-Leu-Ser-Lys-Gly-Cys-Phe-Gly-Leu-Lys-Leu-Asp-Arg-Ile-Gly-Ser- Met-Ser-Gly-Leu-Gly-Cys-OH (Disulfide bridge: 6-22) (3-letter code) GLSKGCFGLKLDRIGSMSGLGC (Disulfide bridge: 6-22) (1-letter code)		
Molecular Weight:	2198.7		
% Peak Area by HPLC	: ≥ 95		
Appearance:	Lyophilized white powder		

Peptide Reconstitution: Using  $H_2O$ , reconstitute by adding 100 µl to 1 mg C-Type Natriuretic peptide.

Storage: C-Type Natriuretic peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at  $-20^{\circ}$ C or lower. Reconstituted peptide can be aliquoted and stored at  $-20^{\circ}$ C or lower.

**Description:** C-type Natriuretic Peptide (CNP), identified in 1990 and called C-type (in order to maintain the alphabetical nomenclature of natriuretic peptides), is the most highly conserved of natriuretic peptides between species. It is derived from a 126 amino acid preprohormone. CNP exists in two mature forms, one found in tissues, another in plasma and cerebrospinal fluid. It is present in high concentration of the vascular tree, especially the endothelium, central nervous tissues, and renal tubular cells. It is a powerful vasorelaxant and inhibitor of smooth muscle cell proliferation and may play a role in coagulation and fibrinogenesis by modulating endothelial cells. Ref: Kalra, P. et al. *Eur. Heart J.* **22**, 997 (2001); Sudoh, T. et al. *Biochem. Biophy. Res. Commun.* **168**, 863 (1990); Koller, KJ. et al. *Science* **252**, 120 (1991).

Additional Information: Listed below are relevant information that may provide a guideline on how to use this product. End users will have to adapt to their own specific applications.

Following a 1.5- to 2-h attachment period, each well was filled with 1 ml of Ham's F-12 culture medium (Gibco-BRL, Gaithersburg, MD); supplemented with 10% fetal bovine serum (Atlanta Biologicals, Atlanta, GA), penicillin (50 U/ml), and streptomycin (50  $\mu$ g/ml); and incubated at 37°C, 5% CO<sub>2</sub>. During the CNP treatment experiments the same culture medium was additionally supplemented with 10<sup>-7</sup> M CNP (32-53; Human, Porcine, AnaSpec, Inc., San Jose, CA). Culture medium was replaced daily-<u>Alan, T. and AC. Tufan *J. Cellular Biochem.* **105**, 227 (2008).</u>

## Published Citations:

Alan, T. and AC. Tufan J. Cellular Biochem. 105, 227 (2008).

## **Related Products:**

Name	<b>Cat #</b>	<b>Size</b>
Atrial Natriuretic Peptide (1-28), human, porcine	AS-20647	0.5 mg
(SLRRSSCFGGRMDRIGAQSGLGCNSFRY (Disulfide bridge: 7-23))	AS-20648	1 mg
Atrial Natriuretic Peptide (1-28), human, porcine, Biotin-labeled (Biotin-SLRRSSCFGGRMDRIGAQSGLGCNSFRY (Disulfide bridge: 7-23))	AS-23972	0.5 mg
Atrial Natriuretic Peptide (1-28), rat	AS-20651	0.5 mg
(SLRRSSCFGGRIDRIGAQSGLGCNSFRY (Disulfide bridge: 7-23))	AS-20652	1 mg
Atrial Natriuretic Peptide (4-18), rat (RSSCFGGRIDRIGAC-NH2 (Disulfide bridge 4-15))	AS-62839	1 mg

For Research Use Only