



Product Data Sheet

Product Name:	Uroguanylin (Rat)	
Catalog Number:	AS-61645 (1mg)	Lot Number: See label on vial
Sequence:	H-Thr-Asp-Glu-Cys-Glu-Leu-Cys-Ile-Asn-Val-Ala-Cys-Thr-Gly-Cys-OH (Disulfide bonds between Cys4-Cys12 and Cys7-Cys15) (3-letter code) TDECELINVACTGC (Disulfide bonds between Cys4-Cys12 and Cys7-Cys15) (1-letter code)	
Molecular Weight:	1569.8	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	

Peptide Reconstitution: Initially try to dissolve Uroguanylin in a small amount of water. If it does not dissolve, use 1% NH₄OH.

Storage: Uroguanylin 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.

Description: Uroguanylin is a natriuretic peptide, a hormone that regulates sodium excretion by the kidney when excess NaCl is consumed. Uroguanylin and guanylin are related peptides that activate common guanylate cyclase signaling molecules in the intestine and kidney. Uroguanylin was isolated from urine and duodenum but was not detected in extracts from the colon of rats. Ref: Fan, X. et al. *Am. J. Physiol. Endocrinol. Metab.* **273**, E957(1997); Forte, L. *J. Clin. Invest.* **112**, 1138 (2003); Jin, X. et al. *Hypertension* **43**, 1133 (2004).

For Research Use Only