



Product Data Sheet

Product Name:	Calcitonin, human	
Catalog Number:	AS-20673 (1 mg) AS-20674 (5 mg)	Lot Number: See label on vial
Sequence:	H-Cys-Gly-Asn-Leu-Ser-Thr-Cys-Met-Leu-Gly-Thr-Tyr-Thr-Gln-Asp-Phe-Asn-Lys-Phe-His-Thr-Phe-Pro-Gln-Thr-Ala-Ile-Gly-Val-Gly-Ala-Pro-NH ₂ (Disulfide bridge: 1-7) (3-letter code) CGNLSTCMLGTYTQDFNKFHTFPQTAIGVGAP-NH ₂ (Disulfide bridge: 1-7) (1-letter code)	
Molecular Weight:	3417.9	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	

Peptide Reconstitution: Calcitonin peptide is freely soluble in water.

Storage: Calcitonin 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: Human calcitonin stimulates cyclic nucleotide accumulation in human kidney cortex and medulla. Calcitonin maybe involved in osteoporosis and in Paget's disease. Ref: Raddino, R. et al. J. Cardiovas. Pharmacol. 29, 463 (1997); Rotella, C. et al. Eur. J. Pharmacol. 107, 347 (1985); Moriarty, et al. Biochem Biophys Res. Commun. 245, 344 (1998); Chen, W. et al. Mol. Pharmacol. 52, 1164 (1997).

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.

Published Citations:

Takiyama, K. et al. *Anal Biochem* **388**, 235 (2009).

For Research Use Only

