

## **Product Data Sheet**

Product Name:	Calcitonin Gene Related Peptide, CGRP, human	
Catalog Number:	AS-20681 (0.5 mg) AS-20682 (1 mg)	Lot Number: See label on vial
Sequence:	ACDTATCVTHRLAGLLSRSGGVVKNNFVPTNVGSKAF-NH2 (Disulfide bridge: 2-7) (3-letter code) ACDTATCVTHRLAGLLSRSGGVVKNNFVPTNVGSKAF-NH2 (Disulfide bridge: 2-7) (1-letter code)	
Molecular Weight:	3789.4	
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: Calcitonin gene-related peptide (CGRP) is a 37-amino acid neuropeptide produced by tissue specific processing of the calcitonin gene and is the major product in neural tissues. CGRP is released by motor neurons where it exerts both short and long term effects on skeletal muscle fibers. In addition, sensory neurons release CGRP to the surrounding vasculature where it is partly responsible for local vasodilation following muscle contraction. CGRP acts through G protein-coupled receptors whose presence and changes in function modulate the peptide<sub>i</sub>'s effects in various tissues. Ref: Nimmagadda D. GS. *Indian J. Pharmacol.* **36**, 277 (2004); Feuerstein, G. et al. *Can. J. Physiol. Pharmacol.* **73**, 1070 (1995).

## Related Products:

Name [Calcitonin Gene Related Peptide, CGRP (8-37), human VTHRLAGLLSRSGGVVKNNFVPTNVGSKAF-NH2	<b>Cat #</b> AS-22857	" <b>Size</b> 1 mg
α-Calcitonin Gene Related Peptide, α-CGRP, rat SCNTATCVTHRLAGLLSRSGGVVKDNFVPTNVGSEAF-NH2	AS-60730	1 mg

For Research Use Only

(Disulfide bridge: 2-7)

©AnaSpec, Inc. 34801 Campus Drive, Fremont, CA 94555 Tel: (800)-452-5530 | service@anaspec.com | www.anaspec.com