



Product Data Sheet

Product Name:	ω-Conotoxin GVIA	
Catalog Number:	AS-22926 (0.5 mg) AS-22927 (1 mg)	Lot Number: See label on vial
Sequence:	H-Cys-Lys-Ser-Hyp-Gly-Ser-Ser-Cys-Ser-Hyp-Thr-Ser-Tyr-Asn-Cys-Cys-Arg-Ser-Cys-Asn-Hyp-Tyr-Thr-Lys-Arg-Cys-Tyr-NH2 (Disulfide bridges: 1-16, 8-19, 15-26) (3-letter code) CKS(Hyp)GSSCS(Hyp)TSYNCCRSCN(Hyp)YTKRCY-NH2 (1-letter code)	
Molecular Weight:	3037.4	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	

Peptide Reconstitution: ω-Conotoxin peptide is freely soluble in water.

Storage: ω-Conotoxin 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: This peptide is a neurotoxin that blocks N-type calcium channels. Ref: Rivier, J. et al. *J. Biol. Chem.* **262**, 1194 (1987); Lampe, RA. et al. *Biochem.* **32**, 3255 (1993); Wright, C. and JA. Angus, *J. Cardiovasc. Pharmacol.* **30**, 392 (1997); Serone, A. and JA. Angus, Br. *J. Pharmacol.* **127**, 927 (1999).

Published Citations:

Baratta, M. V. et al. *J Neurophysiol.* **88**, 3078 (2002).

For Research Use Only