



Product Data Sheet

Product Name:	Charybdotoxin
Catalog Number:	AS-28244 (0.1 mg)
Sequence:	<p>Pyr-Phe-Thr-Asn-Val-Ser-Cys-Thr-Thr-Ser-Lys-Glu-Cys-Trp-Ser-Val-Cys-Gln-Arg-Leu-His-Asn-Thr-Ser-Arg-Gly-Lys-Cys-Met-Asn-Lys-Lys-Cys-Arg-Cys-Tyr-Ser-OH (Disulfide bridge: 7-28, 13-33 and 17-35)</p> <p>Pyr-FTNVSCTTSKECWVQCQLHNTSRGKCMNKKCRCYS (Disulfide bridge: 7-28, 13-33 and 17-35)</p>
Molecular Weight:	4296.1
Peptide Purity:	>95%
Appearance:	Lyophilized white powder
Peptide Reconstitution:	Charybdotoxin peptide is freely soluble in water.
Storage:	Charybdotoxin 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:	Charybdotoxin (ChTX) is a Ca^{2+} -activated K^{+} channel blocker. It depolarizes peripheral T lymphocytes and blocks their mitogen-induced proliferation. ChTX is a highly basic peptide isolated from venom of the scorpion, <i>Leiurus quinquestriatus hebraeus</i> . Ref: Leonard, R. et al. Proc. Natl. Acad. Sci. USA 89, 10094 (1992); Gimenez-Gallego, Proc. Natl. Acad. Sci. USA 85, 3329 (1988). Sugg, E. et al. J. Biol. Chem. 265, 18745 (1990).