



## Product Data Sheet

---

**Product Name:** Chlorotoxin (Cltx)

**Catalog Number:** AS-60770 (0.1 mg)                      **Lot Number:** See label on vial

**Sequence:** H-Met-Cys-Met-Pro-Cys-Phe-Thr-Thr-Asp-His-Gln-Met-Ala-Arg-Lys-Cys-Asp-Asp-Cys-Cys-Gly-Gly-Lys-Gly-Arg-Gly-Lys-Cys-Tyr-Gly-Pro-Gln-Cys-Leu-Cys-Arg-NH<sub>2</sub> (Disulfide bridge: 2-19, 5-28, 16-33, 20-35) (3-letter code)  
MCMPCFTTDHQMARCDDCCGGKGRGKCYGPQCLCR-NH<sub>2</sub> (Disulfide bridge: 2-19, 5-28, 16-33, 20-35) (1-letter code)

**Molecular Weight:** 3996.8

**% Peak Area by HPLC:** ≥ 95

**Appearance:** Lyophilized white powder

**Peptide Reconstitution:** Chlorotoxin peptide is freely soluble in H<sub>2</sub>O.

**Storage:** Chlorotoxin 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:** Chlorotoxin is a 36-amino acid Cl<sup>-</sup> channel blocker from *Leiurus quinquestriatus* scorpion venom. Ref: DeBin, JA. and GR. Strichartz, *Toxicon* **29**, 1403 (1991); DeBin, JA. et al. *Am. J. Physiol.* **264**, C361 (1993).

**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.

Chlorotoxin was obtained from AnaSpec (San Jose, CA, USA). The cells were seeded onto 96-well plates at a density of 6×10<sup>3</sup> cells/well for 24 h (U251-MG and C6 glioma cells) and 7 d (neural cells), respectively. Then SPIO at final concentrations of 5, 10, 25, 50, 75, 100, and 200 mg/L (iron concentrations), chlorotoxin at concentrations of 0.8, 1.6, 2.4, 3.2, 4.0, 4.8, and 5.6 μmol/L, and SPIOFC at final concentrations of 100 mg/L (iron concentrations) and 1.6 μmol/L (chlorotoxin concentration), respectively, were added to the cells and incubated for 24 h-[Meng, X. et al. \*Acta Pharmacol. Sinica\* \*\*28\*\*, 2019 \(2006\).](#)

**Published Citation:**

Meng, X. et al. *Acta Pharmacol. Sinica* **28**, 2019 (2006).

*For Research Use Only*