

PRODUCT DATA SHEET



PEPTIDE NAME: Calciseptine
SEQUENCE: RICYI HKASL PRATK TCVEN TCYKM FIRTQ REYIS ERGCG CPTAM WPYQTECCKG DRCNK

CATALOG NO.: CA-502

LOT NO.: TEMP

MOLECULAR WEIGHT: 7036

PURITY: >98% by HPLC analysis.

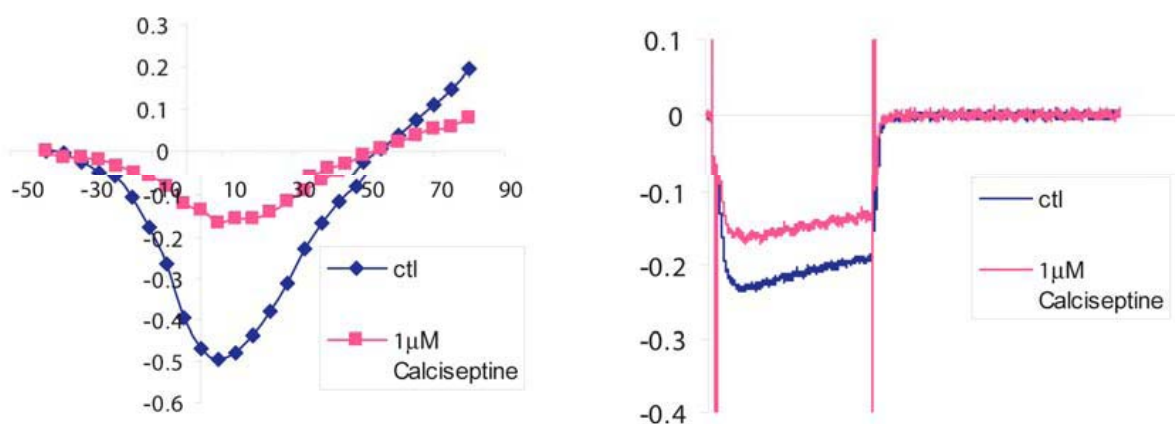
PHYSICAL APPEARANCE: Lyophilized solid.

SOLUBILITY: Any aqueous buffer (pH 7.5).

RECONSTITUTION: Each vial contains 70 µg of unbuffered protein and is provided as a lyophilized powder. Dissolving 70 µg in 10 ml will provide a 1.0 µM stock solution.

Before dissolving the toxin, the tube should be centrifuged. After centrifugation, the toxin must be dissolved into a stock solution using distilled water or an appropriate buffer to a concentration of 10^{-5} - 10^{-6} M. After preparing the stock solution, it should be divided into aliquots and can be stored this way for up to three months at -20°C.

BIOASSAY: Calciseptine was tested for its ability to inhibit L-type Ca^{2+} channels ($Ca_v1.2$, β_3) heterologously expressed in *Xenopus* oocytes. 1 µM Calciseptine blocked 50% of these L-type currents.



Left: I-V relationship. Right: traces of control and inhibited currents.

STORAGE: Lyophilized: 2-3 weeks at room temperature, 1 year at -20°C. Reconstituted: 4 week at 4°C, 3 months at -20°C

REFERENCES:

1. H. Schweitz *et al. Toxicon* 1990 **28** 847
2. J.R. De Weille *et al. Proc. Natl. Acad. Sci (U.S.A.)* 1991 **88** 2437
3. M.C. Garcia *et al J. Membr. Biol.* 2001 **184** 121
4. N. Teramoto *et al. Pflugers Arch.* 1996 **432** 462

HIGHLY TOXIC! USE EXTREME CAUTION! The pharmacological and toxicological properties of this product have not been fully investigated. Avoid all forms of contact with this product by wearing appropriate protective clothing (gloves, mask, lab coat). The Material Safety Data Sheet (MSDS), is available on request. This product must not be used in humans