

PRODUCT DATA SHEET



PEPTIDE NAME: SNX-482
SEQUENCE: GVDKA GCRYM FGGCS VNDDC CPRLG CHSLF SYCAW DLTFD D-

CATALOG NO.: CA-501

LOT NO.: TEMP

MOLECULAR WEIGHT: 4495

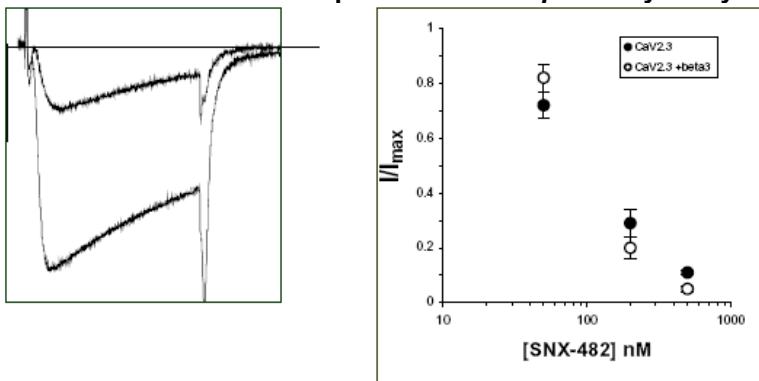
PURITY: >98% by HPLC analysis.

PHYSICAL APPEARANCE: Lyophilized solid.

SOLUBILITY: Any aqueous buffer (pH 7.5).

RECONSTITUTION: Each vial contains 5 µg of unbuffered protein and is provided as a lyophilized powder. Dissolving 5 µg in 1.15 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: Inhibition of Cav2.3 channels expressed in *Xenopus* oocytes by SNX-482.



In *Xenopus* oocytes expressing Cav_v2.3, SNX-482 blocked the currents with an IC₅₀ of about 100 nM.

Left: Block of cloned Cav_v2.3 channel currents (with 10 mM Ba²⁺ as charge carrier) coexpressed with Cav_vβ3 in *Xenopus* oocytes. Currents were elicited by 100 ms pulse from holding potential of -100 mV to 0 mV, every 15 s. Traces show response before and during application of 200 nM SNX-482. The vertical bar represents 1 µA.

Right: Dose response curves for Cav_v2.3 channels with (n=4) or without (n=4) the auxiliary β3 subunit. For these experiments 40 mM Ba²⁺ was used as charge carrier

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. R. Newcomb *et al.* *Biochemistry* 1998 **37** 15353
2. E. Bourmieu *et al.* *Biophys. J.* 2001 **81** 79
3. G. Wang *et al.* *J. Neurosci.* 1999 **19** 9235

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