

PRODUCT INFORMATION



Iberitoxin (trifluoroacetate salt)

Item No. 14608

CAS Registry No.: 182897-30-9

Peptide Sequence: Pyr-FTDVDCSVSKECWSVCKDLFGV
DRGKCMGKKRCRYQ
(Disulfide bridges: 7-28, 13-33, 17-35)

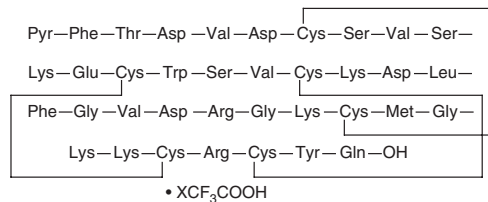
FW: 4,230.9

Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Iberitoxin (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the iberitoxin (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Iberitoxin, a peptide purified from venom of the scorpion *B. tamulus*, is a selective blocker of large-conductance calcium-activated potassium channels ($K_{Ca1.1}/BK_{Ca}$). In single channel recordings, iberitoxin can reversibly block outward currents mediated by $K_{Ca1.1}$ in excised membrane patches from bovine aortic smooth muscle with an IC_{50} value of 250 pM.¹ It is also a noncompetitive, partial inhibitor of charybdotoxin binding in bovine aortic sarcolemmal membrane vesicles ($K_i = 250$ pM).¹

Reference

- Galvez, A., Gimenez-Gallego, G., Reuben, J.P., *et al.* Purification and characterization of a unique, potent, peptidyl probe for the high conductance calcium-activated potassium channel from venom of the scorpion *Buthus tamulus*. *J. Biol. Chem.* **265**(19), 11083-11090 (1990).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/02/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM