

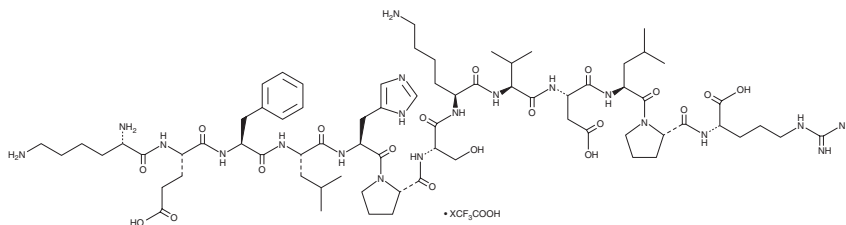
PRODUCT INFORMATION



SOR-C13 (trifluoroacetate salt)

Item No. 36764

Formal Name: L-lysyl-L- α -glutamyl-L-phenylalanyl-L-leucyl-L-histidyl-L-prolyl-L-seryl-L-lysyl-L-valyl-L- α -aspartyl-L-leucyl-L-prolyl-L-arginine, trifluoroacetate salt
MF: C₇₂H₁₁₆N₂₀O₁₉ • XCF₃COOH
FW: 1,565.8
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

SOR-C13 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the SOR-C13 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. SOR-C13 (trifluoroacetate salt) is soluble in the organic solvent DMSO at a concentration of approximately 12 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of SOR-C13 (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of SOR-C13 (trifluoroacetate salt) in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

SOR-C13 is a peptide antagonist of transient receptor potential vanilloid 6 (TRPV6).¹ It inhibits calcium flux in HEK293 cells expressing human TRPV6 (EC₅₀ = 14 nM). SOR-C13 (400, 600, and 800 mg/kg) reduces tumor volume in an SKOV3 ovarian cancer mouse xenograft model.

Reference

1. Xue, H., Wang, Y., MacCormack, T.J., *et al.* Inhibition of Transient Receptor Potential Vanilloid 6 channel, elevated in human ovarian cancers, reduces tumour growth in a xenograft model. *J. Cancer* **9**(17), 3196-3207 (2018).

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, 12/21/2022

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