

PRODUCT INFORMATION



Caloxin 2A1 (trifluoroacetate salt)

Item No. 29956

Formal Name: L-valyl-L-seryl-L-asparaginyl-L-seryl-L-asparaginyl-L-tryptophyl-L-prolyl-L-seryl-L-phenylalanyl-L-prolyl-L-seryl-L-serylglycylglycyl-glycinamide, trifluoroacetate salt

Synonym: VSNSNWPSFPSSGGG-NH₂

MF: C₆₄H₉₁N₁₉O₂₂ • XCF₃COOH

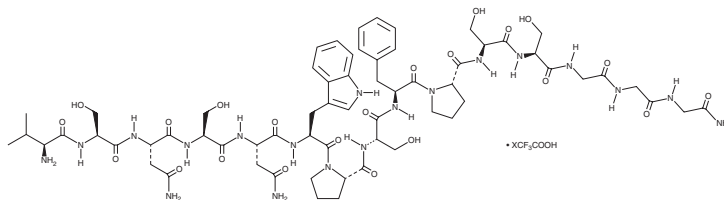
FW: 1,478.5

Purity: ≥98%

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

Caloxin 2A1 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the caloxin 2A1 (trifluoroacetate salt) in water. The solubility of caloxin 2A1 (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Caloxin 2A1 is a peptide plasma membrane Ca²⁺/Mg²⁺-ATPase inhibitor (IC₅₀ = 0.4 mM in erythrocyte leaky ghosts).¹ It is selective for the plasma membrane Ca²⁺/Mg²⁺-ATPase over the erythrocyte ghost Mg²⁺- and Na⁺/K⁺-ATPases and the skeletal muscle sarcoplasmic reticulum Ca²⁺/Mg²⁺-ATPase at 0.9 mM. Caloxin 2A1 (0.34 mM) induces relaxation in precontracted rat aortic rings, an effect that can be reversed by the nitric oxide synthase inhibitor L-NAME (Item No. 80210).

Reference

1. Chaudhary, J., Walia, M., Matharu, J., *et al.* Caloxin: A novel plasma membrane Ca²⁺ pump inhibitor. *Am. J. Physiol. Cell Physiol.* **280**(4), C1027-C1030 (2001).

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, 10/31/2022

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