

PRODUCT INFORMATION



Calcineurin Substrate (trifluoroacetate salt)

Item No. 34670

Formal Name: L- α -aspartyl-L-leucyl-L- α -aspartyl-L-valyl-L-prolyl-L-isoleucyl-L-prolylglycyl-L-arginyl-L-phenylalanyl-L- α -aspartyl-L-arginyl-L-arginyl-L-valyl-L-seryl-L-valyl-L-alanyl-L-alanyl-L-glutamic acid, trifluoroacetate salt

Synonyms: H-Asp-Leu-Asp-Val-Pro-Ile-Pro-Gly-Arg-Phe-Asp-Arg-Arg-Val-Ser-Val-Ala-Ala-Glu-OH, H-DLDVPIPGRFDRRVSVAEE-OH

MF: C₉₂H₁₅₀N₂₈O₂₉ • XCF₃COOH

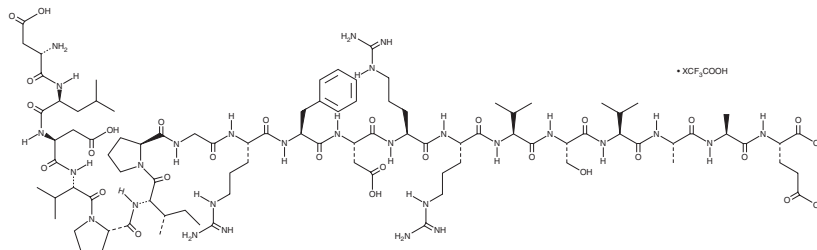
FW: 2,112.4

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

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

Description

Calcineurin substrate is a synthetic peptide that corresponds to the regulatory RII subunit of protein kinase A (PKA).¹ It can be phosphorylated by PKA, and phosphorylated calcineurin substrate can be used to quantify calcineurin activity.

Reference

1. Namgaladze, D., Hofer, H.W., and Ullrich, V. Redox control of calcineurin by targeting the binuclear Fe²⁺-Zn²⁺ center at the enzyme active site. *J. Biol. Chem.* **277**(8), 5962-5969 (2002).

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, 11/22/2022

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