

PRODUCT INFORMATION



DAPK Substrate Peptide (trifluoroacetate salt)

Item No. 35221

Formal Name: L-lysyl-L-lysyl-L-arginyl-L-prolyl-L-glutaminy-L-arginyl-L-arginyl-L-tyrosyl-L-seryl-L-asparaginy-L-valyl-L-phenylalanine, trifluoroacetate Substrate Peptide

Synonym: Death-associated Protein Kinase Substrate Peptide

MF: C₇₀H₁₁₅N₂₅O₁₇ • XCF₃COOH

FW: 1,578.8

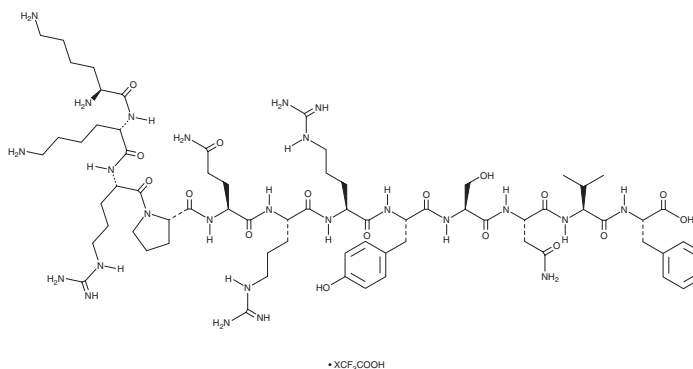
Purity: ≥98%

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

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

Description

DAPK substrate peptide is a peptide substrate for death-associated protein kinase 1 (DAPK1), a serine/threonine kinase involved in apoptosis and tumor suppression.¹ It has been used as a target substrate for the quantification of DAPK1 activity.

Reference

1. Velentza, A.V., Schumacher, A.M., Weiss, C., *et al.* A protein kinase associated with apoptosis and tumor suppression: Structure, activity, and discovery of peptide substrates. *J. Biol. Chem.* **276(42)**, 38956-38965 (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, 12/12/2022

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