

PRODUCT INFORMATION



R18 (trifluoroacetate salt)

Item No. 36695

Formal Name: L-prolyl-L-histidyl-L-cysteinyl-L-valyl-L-prolyl-L-arginyl-L- α -aspartyl-L-leucyl-L-seryl-L-tryptophyl-L-leucyl-L- α -aspartyl-L-leucyl-L- α -glutamyl-L-alanyl-L-asparaginyl-L-methionyl-L-cysteinyl-L-leucyl-L-proline, trifluoroacetate salt

Peptide Sequence: PHCVPRDLSWLDLEANMCLP-OH

MF: C₁₀₁H₁₅₇N₂₇O₂₉S₃ • XCF₃COOH

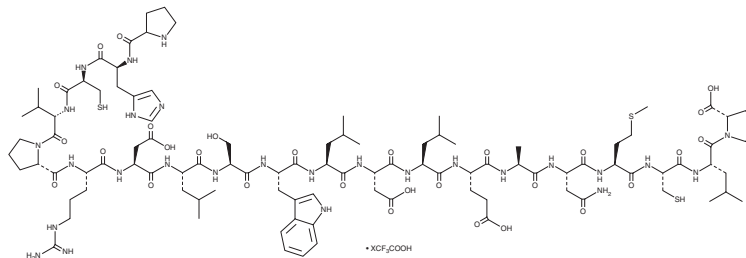
FW: 2,309.7

Purity: \geq 95%

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

R18 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the R18 (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

R18 is a synthetic peptide inhibitor of the protein-protein interaction between 14-3-3 proteins and Raf-1.¹ It binds to 14-3-3 τ , - β , and - ζ (K_d s = 7.1, 7, and 8.9 nM, respectively) and inhibits 14-3-3 τ binding to Raf-1 in a filter binding assay when used at a concentration of 5 nM. R18 also inhibits 14-3-3 ζ -mediated protection of Raf-1 from phosphatase-induced inactivation.

Reference

1. Wang B, Yang H, Liu Y.C., *et al.* Isolation of high-affinity peptide antagonists of 14-3-3 proteins by phage display. *Biochemistry* **38**(38), 12499-12504 (1999).

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, 07/24/2023

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