

PRODUCT INFORMATION

RS Repeat Peptide (trifluoroacetate salt)

Item No. 37498

Formal Name: glycyl-L-arginyl-L-seryl-L-arginyl-L-seryl-L-arginyl-L-seryl-L-arginyl-L-seryl-L-arginyl-L-seryl-L-arginine, trifluoroacetate salt

Peptide Sequence: GRSRSRSRSRSRSR-OH

MF: $C_{71}H_{136}N_{40}O_{24} \bullet XCF_3COOH$

FW: 1,934.1

Purity:

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

H-Gly-Arg-Ser-Arg-Ser-Arg-Ser-Arg-Ser-Arg-

$$\text{Ser}-\text{Arg}-\text{Ser}-\text{Arg}-\text{Ser}-\text{Arg}-\text{OH}$$

- XCF_3COOH

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

Laboratory Procedures

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

Description

RS repeat peptide is a serine/arginine-rich peptide substrate for CDC-like kinase 1 (Clk1).^{1,2} It has been used to identify inhibitors of Clk1 with anticancer activity.

References

1. Schmitt, C., Kail, D., Mariano, M., *et al.* Design and synthesis of a library of lead-like 2,4-bisheterocyclic substituted thiophenes as selective Dyrk/Clk inhibitors. *PLoS One* **9**(3), e87851 (2014).
2. Araki, S., Dairiki, R., Nakayama, Y., *et al.* Inhibitors of CLK protein kinases suppress cell growth and induce apoptosis by modulating pre-mRNA splicing. *PLoS One* **10**(1), e0116929 (2015).

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, 01/09/2024

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