

PRODUCT INFORMATION

Histone H1-derived Peptide (trifluoroacetate salt)

Item No. 27439

Formal Name: glycylglycylglycyl-L-prolyl-L-alanyl-L-threonyl-L-prolyl-L-lysyl-L-lysyl-L-alanyl-L-lysyl-L-lysyl-L-leucine, trifluoroacetate salt

Synonym: GGGPATPKKAKKL H—Gly—Gly—Gly—Pro—Ala—Thr—Pro—Lys—Lys—Ala—Lys—Lys—Leu—OH

MF: $C_{56}H_{101}N_{17}O_{15} \cdot XCF_3COOH$

FW: 1,252.5

Purity: $\geq 95\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years

$\bullet XCF_3COOH$

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

Laboratory Procedures

Histone H1-derived peptide (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the histone H1-derived peptide (trifluoroacetate salt) in water. The solubility of histone H1-derived peptide (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Histone H1-derived peptide is a substrate for the *Leishmania* complex between Cdc2-related kinase 3 (CRK3) and cyclin CYC6 (CRK3/CYC6).¹ It is phosphorylated by CRK3/CYC6 and has been used in high-throughput screening assays for the identification of CRK3/CYC6 inhibitors.

Reference

1. Walker, R.G., Thomson, G., Malone, K., et al. High throughput screens yield small molecule inhibitors of *Leishmania* CRK3:CYC6 cyclin-dependent kinase. *PLoS Negl. Trop. Dis.* **5**(4), e1033 (2011).

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, 11/30/2022

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