

PRODUCT INFORMATION



Cdk2 Peptide Substrate (trifluoroacetate salt)

Item No. 44656

Formal Name: L-histidyl-L-histidyl-L-alanyl-L-seryl-L-prolyl-L-arginyl-L-lysine, trifluoroacetate salt

Synonyms: Cyclin-dependent Kinase 2 Peptide Substrate, His-His-Ala-Ser-Pro-Arg-Lys-OH

Peptide Sequence: HHASPRK-OH

MF: C₃₅H₅₇N₁₅O₉ • XCF₃COOH

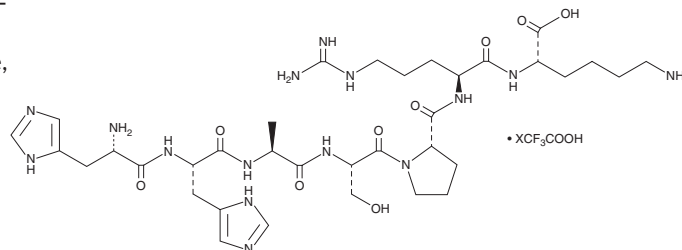
FW: 831.9

Purity: ≥90%

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

Cdk2 peptide substrate (trifluoroacetate salt) is supplied as a solid. Aqueous solutions of Cdk2 peptide substrate (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. Cdk2 peptide substrate (trifluoroacetate salt) is soluble (≥10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Cdk2 peptide substrate is an optimal peptide substrate of Cdk2.¹ It binds across the Cdk2 catalytic site when in complex with Cdk2 phosphorylated at T160 and cyclin A3. Cdk2 peptide substrate has been used in enzyme assays for Cdk2 activity.^{2,3}

References

1. Brown, N.R., Noble, M.E., Endicott, J.A., *et al.* The structural basis for specificity of substrate and recruitment peptides for cyclin-dependent kinases. *Nat. Cell. Biol.* **1(7)**, 438-443 (1999).
2. Anscombe, E., Meschini, E., Mora-Vidal, R., *et al.* Identification and characterization of an irreversible inhibitor of CDK2. *Chem. Biol.* **22(9)**, 1159-1164 (2015).
3. Song, K., Liu, X., Huang, W., *et al.* Improved method for the identification and validation of allosteric sites. *J. Chem. Inf. Model* **57(9)**, 2358-2363 (2017).

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, 04/14/2026

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