

PRODUCT INFORMATION



Dnp-PLG-Cys(Me)-HA-DArg-NH₂ (trifluoroacetate salt)

Item No. 24663

CAS Registry No.: 145224-98-2
Formal Name: 1-(2,4-dinitrophenyl)-L-prolyl-L-leucylglycyl-S-methyl-L-cysteinyl-L-histidyl-L-alanyl-D-argininamide, trifluoroacetate salt

Synonyms: Dnp-Pro-Leu-Gly-Cys(Me)-His-Ala-DArg-NH₂, Matrix Metalloproteinase-1 Substrate II, MMP-1 Substrate II

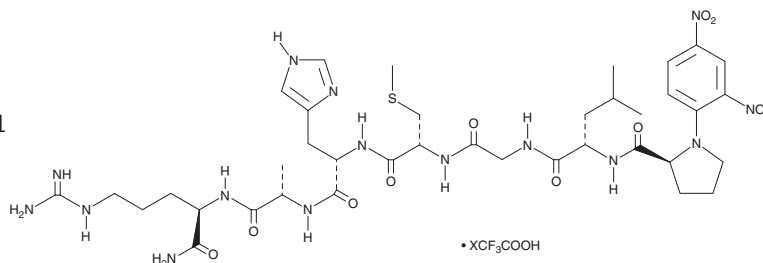
MF: C₃₈H₅₇N₁₅O₁₁S • XCF₃COOH
FW: 932.0

Purity: ≥95%

Supplied as: A lyophilized powder

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

Dnp-PLG-Cys(Me)-HA-DArg-NH₂ (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the Dnp-PLG-Cys(Me)-HA-DArg-NH₂ (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Dnp-PLG-Cys(Me)-HA-DArg-NH₂ (trifluoroacetate salt) is soluble in the organic solvent formic acid at a concentration of approximately 1 mg/ml.

Description

Dnp-PLG-Cys(Me)-HA-DArg-NH₂ is a peptide substrate for matrix metalloproteinase-1 (MMP-1).¹ It has an improved k_{cat}/K_m ratio compared with the MMP-1 fluorogenic substrate Dnp-PLGLWA-DArg-NH₂.

Reference

1. Berman, J., Green, M., Sugg, E., *et al.* Rapid optimization of enzyme substrates using defined substrate mixtures. *J. Biol. Chem.* **267**(3), 1434-1437 (1992).

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, 12/17/2022

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