

PRODUCT INFORMATION



Dnp-PLGMWSR (trifluoroacetate salt)

Item No. 24698

Formal Name: 1-(2,4-dinitrophenyl)-L-prolyl-L-leucylglycyl-L-methionyl-L-tryptophyl-L-seryl-L-arginine, 2,2,2-trifluoroacetate

Synonyms: Dnp-Pro-Leu-Gly-Met-Trp-Ser-Arg-OH, Matrix Metalloproteinase-2/Matrix Metalloproteinase-9 Fluorogenic Substrate I, MMP-2/MMP-9 Fluorogenic Substrate I

MF: C₄₄H₆₁N₁₃O₁₃S • XCF₃COOH

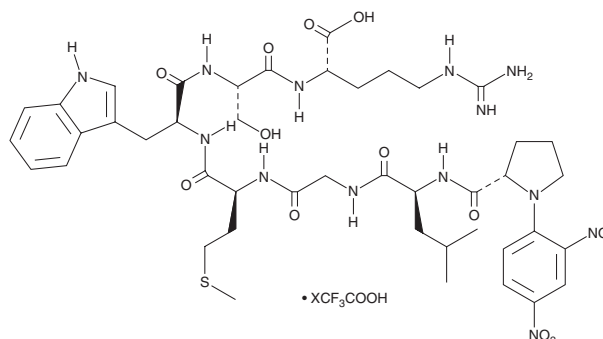
FW: 1,012.1

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

Description

Dnp-PLGMWSR is a fluorogenic substrate for matrix metalloproteinase-2 (MMP-2) and MMP-9.¹ The activity of MMP-2 and MMP-9 can be quantified by measuring tryptophan fluorescence that is unquenched upon peptide hydrolysis that removes the N-terminal dinitrophenol (Dnp) group.

Reference

1. Netzel-Arnett, S., Mallya, S.K., Nagase, H., *et al.* Continuously recording fluorescent assays optimized for five human matrix metalloproteinases. *Anal. Biochem.* **195**(1), 86-92 (1991).

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, 12/01/2022

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