

PRODUCT INFORMATION



Dnp-GPLGMRGL-NH₂ (trifluoroacetate salt)

Item No. 24701

Formal Name: N-(2,4-dinitrophenyl)glycyl-L-prolyl-L-leucylglycyl-L-methionyl-L-arginylglycyl-L-leucinamide, 2,2,2-trifluoroacetate

Synonyms: Dnp-Gly-Pro-Leu-Gly-Met-Arg-Gly-Leu-NH₂, Matrix Metalloproteinase-13 Substrate, MMP-13 Substrate

MF: C₄₀H₆₄N₁₄O₁₂S • XCF₃COOH

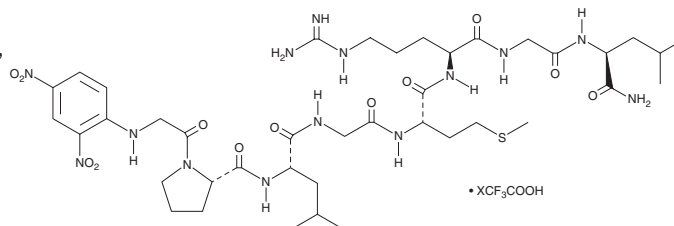
FW: 965.1

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

Description

Dnp-GPLGMRGL-NH₂ is a peptide substrate for matrix metalloproteinase-13 (MMP-13) with a k_{cat}/K_m value of 4,200,000 M⁻¹s⁻¹.¹

Reference

1. Deng, S.-J., Bickett, D.M., Mitchell, J.L., *et al.* Substrate specificity of human collagenase 3 assessed using a phage-displayed peptide library. *J. Biol. Chem.* **275**(40), 31422-31427 (2000).

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, 11/17/2022

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