

PRODUCT INFORMATION



Dnp-GPLG

Item No. 24705

Formal Name: (2,4-dinitrophenyl)glycyl-L-prolyl-L-leucylglycine
Synonyms: Dnp-Gly-Pro-Leu-Gly-OH,
Matrix Metalloproteinase Substrate II Control,
MMP Substrate II Control

MF: C₂₁H₂₈N₆O₉

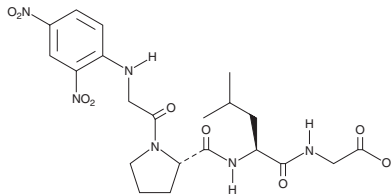
FW: 508.5

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

Description

Dnp-GPLG is a peptide for use as a negative control for activity of matrix metalloproteinases (MMPs) that cleave peptides containing the sequence Dnp-GPLG.

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