

# PRODUCT INFORMATION

## Dnp-P-Cha-G-Cys(Me)-HA-K(Nma)-NH<sub>2</sub> (trifluoroacetate salt)

Item No. 24709

**Formal Name:** 1-(2,4-dinitrophenyl)-L-prolyl-3-cyclohexyl-L-alanylglycyl-S-methyl-L-cysteiny-L-histidyl-L-alanyl-N<sup>6</sup>-[2-(methylamino)benzoyl]-L-lysineamide, trifluoroacetate salt

**Synonyms:** Dnp-Pro-Cha-Gly-Cys(Me)-His-Ala-Lys(Nma)-NH<sub>2</sub>, Matrix Metalloproteinase-1/Matrix Metalloproteinase-9 Fluorogenic Substrate, MMP-1/MMP-9 Fluorogenic Substrate

**MF:** C<sub>49</sub>H<sub>68</sub>N<sub>14</sub>O<sub>12</sub>S • XCF<sub>3</sub>COOH

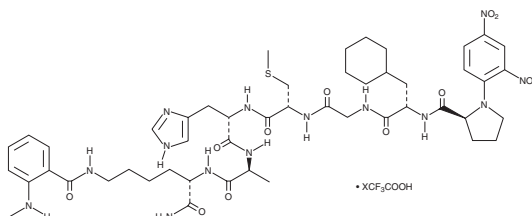
**FW:** 1,077.2

**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-P-Cha-G-Cys(Me)-HA-K(Nma)-NH<sub>2</sub> (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the Dnp-P-Cha-G-Cys(Me)-HA-K(Nma)-NH<sub>2</sub> (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Dnp-P-Cha-G-Cys(Me)-HA-K(Nma)-NH<sub>2</sub> (trifluoroacetate salt) is soluble in the organic solvent formic acid at a concentration of approximately 1 mg/ml.

### Description

Dnp-P-Cha-G-Cys(Me)-HA-K(Nma)-NH<sub>2</sub> is a fluorogenic substrate for matrix metalloproteinase-1 (MMP-1) and MMP-9.<sup>1</sup> Upon cleavage by MMP-1 or MMP-9, N-methylantranilic acid (Nma) is unquenched and its fluorescence can be used to quantify MMP activity. Nma displays excitation/emission spectra of 340/440 nm, respectively.

### Reference

1. Bickett, D.M., Green, M.D., Berman, J., *et al.* A high throughput fluorogenic substrate for interstitial collagenase (MMP-1) and gelatinase (MMP-9). *Anal. Biochem.* **212(1)**, 58-64 (1993).

#### 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, 08/17/2023

#### 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