

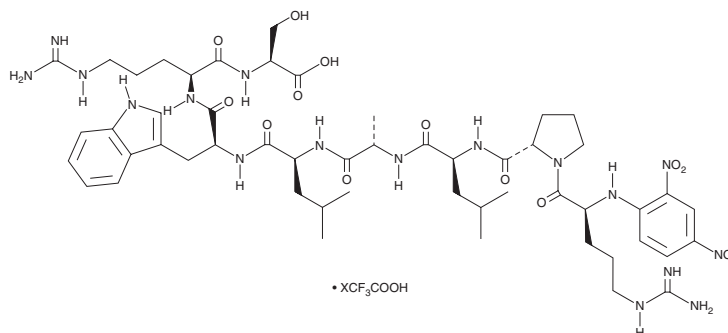
# PRODUCT INFORMATION



## Dnp-RPLALWRS (trifluoroacetate salt)

Item No. 24699

**Formal Name:** N<sup>2</sup>-(2,4-dinitrophenyl)-L-arginyl-L-prolyl-L-leucyl-L-alanyl-L-leucyl-L-tryptophyl-L-arginyl-L-serine, 2,2,2-trifluoroacetate  
**Synonyms:** Dnp-Arg-Pro-Leu-Ala-Leu-Trp-Arg-Ser-OH, Matrix Metalloproteinase-7 Fluorogenic Substrate, MMP-7 Fluorogenic Substrate  
**MF:** C<sub>52</sub>H<sub>77</sub>N<sub>17</sub>O<sub>14</sub> • XCF<sub>3</sub>COOH  
**FW:** 1,164.3  
**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-RPLALWRS (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the Dnp-RPLALWRS (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Dnp-RPLALWRS (trifluoroacetate salt) is soluble in the organic solvent formic acid at a concentration of approximately 1 mg/ml.

### Description

Dnp-RPLALWRS is a fluorogenic substrate for matrix metalloproteinase-7 (MMP-7).<sup>1</sup> The activity of MMP-7 can be quantified by measuring tryptophan fluorescence that is unquenched upon peptide hydrolysis that removes the N-terminal dinitrophenol (Dnp) group.

### Reference

1. Welch, A.R., Holman, C.N., Browner, M.F., *et al.* Purification of human matrilysin produced in *Escherichia coli* and characterization using a new optimized fluorogenic peptide substrate. *Arch. Biochem. Biophys.* **324**(1), 59-64 (1995).

#### 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