

# PRODUCT INFORMATION



## Abz-Val-Ala-Asp-Nva-Arg-Asp-Arg-Gln-EDDnp (trifluoroacetate salt)

Item No. 9002021

**Formal Name:** N-(2-aminobenzoyl)-L-valyl-L-alanyl-L- $\alpha$ -aspartyl-L-norvalyl-L-arginyl-L- $\alpha$ -aspartyl-L-arginyl-N<sup>1</sup>-[2-[(2,4-dinitrophenyl)amino]ethyl]-L-glutamamide, trifluoroacetate salt

**MF:** C<sub>53</sub>H<sub>80</sub>N<sub>20</sub>O<sub>18</sub> • XCF<sub>3</sub>COOH

**FW:** 1,285.3

**Purity:** ≥98%

**UV/Vis.:**  $\lambda_{\max}$ : 252, 345 nm

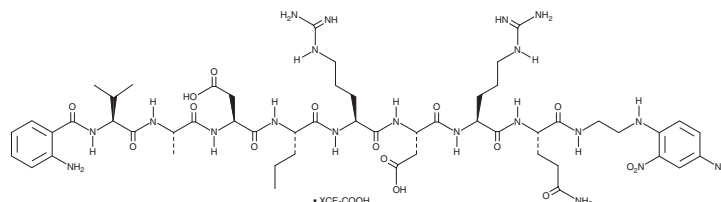
**Supplied as:** A lyophilized powder

**Storage:** -20°C

**Stability:** ≥4 years

**Special Conditions:** Light sensitive

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.



### Laboratory Procedures

Abz-Val-Ala-Asp-Nva-Arg-Asp-Arg-Gln-EDDnp (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the Abz-Val-Ala-Asp-Nva-Arg-Asp-Arg-Gln-EDDnp (trifluoroacetate salt) in the solvent of choice. Abz-Val-Ala-Asp-Nva-Arg-Asp-Arg-Gln-EDDnp (trifluoroacetate salt) is soluble in organic solvents such as DMSO which should be purged with an inert gas. Abz-Val-Ala-Asp-Nva-Arg-Asp-Arg-Gln-EDDnp (trifluoroacetate salt) is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

### Description

Abz-Val-Ala-Asp-Nva-Arg-Asp-Arg-Gln-EDDnp is a fluorescence-quenched peptide substrate for human proteinase 3 ( $k_{\text{cat}}/K_m = 1,570 \text{ mM}^{-1}\text{s}^{-1}$ ).<sup>1</sup> Enzymatic peptide hydrolysis disrupts the Abz:EDDnp donor-acceptor pair allowing fluorescence measurement of Abz (ex = 320 nm; em = 420 nm).

### Reference

1. Kormaz, B., Attucci, S., Juliano, M.A., *et al.* Measuring elastase, proteinase 3 and cathepsin G activities at the surface of human neutrophils with fluorescence resonance energy transfer substrates. *Nat. Protoc.* **3**(6), 991-1000 (2008).

#### 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, 12/13/2022

#### 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