

# PRODUCT INFORMATION



## Z-VDVAD-pNA (trifluoroacetate salt)

Item No. 27426

**Formal Name:** N-[(phenylmethoxy)carbonyl]-L-valyl-L- $\alpha$ -aspartyl-L-valyl-L-alanyl-N-(4-nitrophenyl)-L- $\alpha$ -asparagine, trifluoroacetate salt

**Synonyms:** Z-VDVAD-*p*-nitroanilide,  
Z-Val-Asp-Val-Ala-Asp-*p*NA,  
Z-Val-Asp-Val-Ala-Asp-*p*-nitroanilide

**MF:** C<sub>35</sub>H<sub>45</sub>N<sub>7</sub>O<sub>13</sub> • XCF<sub>3</sub>COOH

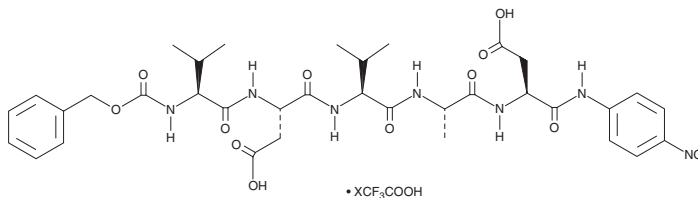
**FW:** 771.8

**Purity:** ≥95%

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥4 years



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

### Laboratory Procedures

Z-VDVAD-*p*NA (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the Z-VDVAD-*p*NA (trifluoroacetate salt) in the solvent of choice. Z-VDVAD-*p*NA (trifluoroacetate salt) is soluble in 5% NH<sub>3</sub> (aqueous). The solubility of Z-VDVAD-*p*NA (trifluoroacetate salt) in 5% NH<sub>3</sub> (aqueous) is approximately 1 mg/ml.

### Description

Z-VDVAD-*p*NA is a colorimetric substrate for caspase-2.<sup>1</sup> Caspase-2 binds to and cleaves the Val-Asp-Val-Ala-Asp (VDVAD) peptide sequence to release *p*-nitroanilide (*p*NA), which can be quantified by colorimetric detection at 405 nm as a measure of caspase-2 activity. Z-VDVAD-*p*NA can also be cleaved, with lower efficiency, by caspase-3 and -7.

### Reference

1. Talanian, R.V., Quinlan, C., Trautz, S., *et al.* Substrate specificities of caspase family proteases. *J. Biol. Chem.* **272**(15), 9677-9682 (1997).

#### 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, 12/08/2022

#### 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](http://WWW.CAYMANCHEM.COM)