

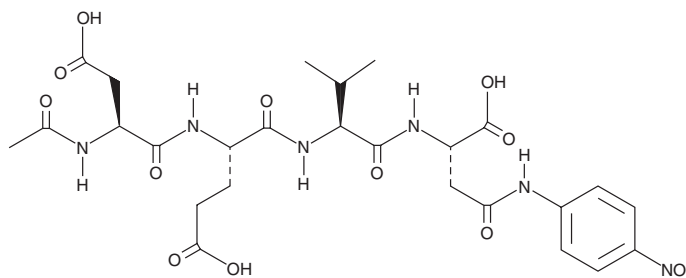
# PRODUCT INFORMATION



## Ac-DEVD-pNA

Item No. 14460

**CAS Registry No.:** 189684-50-2  
**Formal Name:** N-acetyl-L- $\alpha$ -aspartyl-L- $\alpha$ -glutamyl-L-valyl-N-(4-nitrophenyl)-L- $\alpha$ -asparagine  
**Synonym:** Ac-Asp-Glu-Val-Asp-pNA  
**MF:** C<sub>26</sub>H<sub>34</sub>N<sub>6</sub>O<sub>13</sub>  
**FW:** 638.6  
**Purity:**  $\geq$ 95%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years



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

### Laboratory Procedures

Ac-DEVD-pNA is supplied as a crystalline solid. Ac-DEVD-pNA is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, Ac-DEVD-pNA should first be dissolved in methanol and then diluted with the aqueous buffer of choice. Ac-DEVD-pNA has a solubility of approximately 1 mg/ml in a 1:1 solution of acetonitrile:methanol (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

During apoptosis, activated caspase-3 cleaves several substrates, including poly(ADP-ribose) polymerase, which it specifically targets at the amino sequence DEVD.<sup>1</sup> Ac-DEVD-pNA is a *para*-nitro aniline chromophore cleaved by caspases ( $K_m$  = 18, 11, 32, 180, and 12  $\mu$ M for caspases-1, -3, -4, -6, and -7, respectively).<sup>2</sup> It is not cleaved by caspase-2.<sup>2</sup> Cleavage is monitored colorimetrically at 405 nm.

### References

1. Lazebnik, Y.A., Kaufmann, S.H., Desnoyers, S., *et al.* Cleavage of poly(ADP-ribose) polymerase by a proteinase with properties like ICE. *Nature* **371**, 346-347 (1994).
2. Talanian, R.V., Quinlan, C., Trautz, S., *et al.* Substrate specificities of caspase family proteases. *J. Biol. Chem.* **272**, 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, 11/04/2022

#### 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