

PRODUCT INFORMATION



Ac-VEID-pNA (trifluoroacetate salt)

Item No. 27104

Formal Name: N-acetyl-L-valyl-L- α -glutamyl-L-isoleucyl-N-(4-nitrophenyl)-L- α -asparagine, trifluoroacetate salt

Synonyms: Ac-Val-Glu-Ile-Asp-pNA, Ac-Val-Glu-Ile-Asp-p-nitroanilide, Caspase-6 Chromogenic Substrate

MF: C₂₈H₄₀N₆O₁₁ • XCF₃COOH

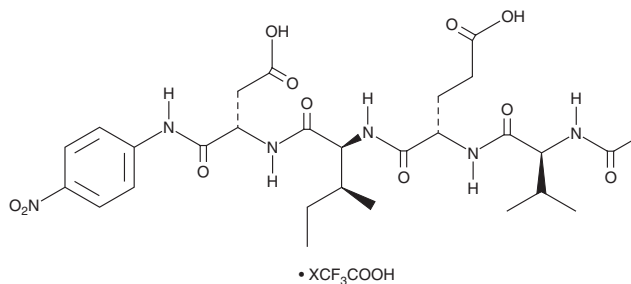
FW: 636.7

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

Ac-VEID-pNA (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the Ac-VEID-pNA (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Ac-VEID-pNA (trifluoroacetate salt) is soluble in the organic solvent DMSO. The solubility of Ac-VEID-pNA (trifluoroacetate salt) in DMSO is approximately 1 mg/ml.

Aqueous solutions of Ac-VEID-pNA (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of Ac-VEID-pNA (trifluoroacetate salt) in 5% NH₃ is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Ac-VEID-pNA is a substrate for caspase-6. Caspase-6 binds to and cleaves the Val-Glu-Ile-Asp (VEID) peptide sequence to release p-nitroanilide (pNA), which can be quantified by colorimetric detection at 405 nm as a measure of enzyme activity.¹

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, 12/05/2022

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