

PRODUCT INFORMATION



Ac-IEPD-*p*NA (trifluoroacetate salt)

Item No. 27139

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

Synonyms: Ac-Ile-Glu-Pro-Asp-*p*-nitroanilide, Caspase-8 Chromogenic Substrate, Granzyme B Substrate VIII

MF: C₂₈H₃₈N₆O₁₁ • XCF₃COOH

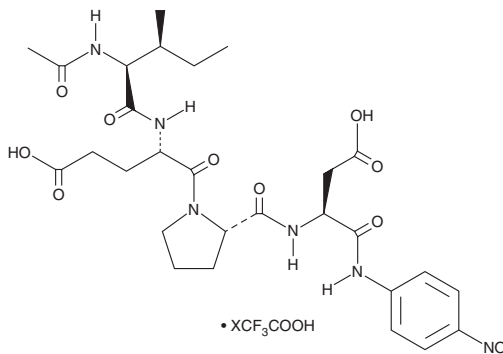
FW: 634.6

Purity: \geq 95%

Supplied as: A 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-IEPD-*p*NA (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the Ac-IEPD-*p*NA (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Ac-IEPD-*p*NA (trifluoroacetate salt) is soluble in the organic solvent DMSO at a concentration of approximately 1 mg/ml.

Description

Ac-IEPD-*p*NA is a colorimetric substrate for granzyme B and caspase-8.^{1,2} Granzyme B and caspase-8 preferentially bind to and cleave the Ile-Glu-Pro-Asp (IEPD) peptide sequence to release *p*-nitroanilide (*p*NA), which can be quantified by colorimetric detection at 405 nm as a measure of granzyme B and caspase-8 activity.

References

1. Ewen, C., Kane, K.P., Shostak, I., *et al.* A novel cytotoxicity assay to evaluate antigen-specific CTL responses using a colorimetric substrate for Granzyme B. *J. Immunol. Methods* **276**(1-2), 89-101 (2003).
2. Almeida, S., Domingues, A., Rodrigues, L., *et al.* FK506 prevents mitochondrial-dependent apoptotic cell death induced by 3-nitropropionic acid in rat primary cortical cultures. *Neurobiol. Dis.* **17**(3), 435-444 (2004).

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, 11/07/2022

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