

PRODUCT INFORMATION

Ac-IETD-CHO (trifluoroacetate salt)

Item No. 27100

Formal Name: N-acetyl-L-isoleucyl-L- α -glutamyl-N-[(1S)-2-carboxy-1-formylethyl]-L-threoninamide, trifluoroacetate salt

Synonyms: Ac-Ile-Glu-Thr-Asp-CHO, Caspase-8 Inhibitor

MF: C₂₁H₃₄N₄O₁₀ • XCF₃COOH

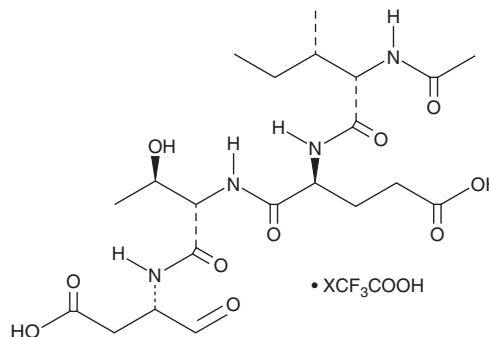
FW: 502.5

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-IETD-CHO (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the Ac-IETD-CHO (trifluoroacetate salt) in water. The solubility of Ac-IETD-CHO (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Ac-IETD-CHO is an inhibitor of caspase-8 (IC₅₀ = 5 nM).¹ It inhibits fluticasone propionate-induced increases in apoptosis in peripheral blood T-cells isolated from patients with asthma when used at a concentration 60 μ M.² Ac-IETD-CHO (10 mg/kg) inhibits increases in hepatic caspase-3 activity, DNA fragmentation, cytosolic cytochrome c levels, and proteolytic processing of caspase-3, -8, and -9 proenzymes in a mouse model of Fas-induced apoptosis and liver failure.³

References

1. Smith, G.K., Barrett, D.G., Blackburn, K.J., *et al.* Expression, preparation, and high-throughput screening of caspase-8: Discovery of redox-based and steroid diacid inhibition. *Arch. Biochem. Biophys.* **399**(2), 195-205 (2002).
2. Pace, E., Gagliardo, R., Melis, M., *et al.* Synergistic effects of fluticasone propionate and salmeterol on *in vitro* T-cell activation and apoptosis in asthma. *J. Allergy. Clin. Immunol.* **114**(5), 1216-1223 (2004).
3. Bajt, M.L., Lawson, J.A., Vonderfecht, S.L., *et al.* Protection against Fas receptor-mediated apoptosis in hepatocytes and nonparenchymal cells by a caspase-8 inhibitor *in vivo*: Evidence for a postmitochondrial processing of caspase-8. *Toxicol. Sci.* **58**(1), 109-117 (2000).

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, 11/11/2022

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