

PRODUCT INFORMATION



Ac-WEAD-pNA (trifluoroacetate salt)

Item No. 27424

Formal Name: (4S,7S,10S,13S)-4-((1H-indol-3-yl)methyl)-7-(2-carboxyethyl)-10-methyl-13-((4-nitrophenyl)carbamoyl)-2,5,8,11-tetraoxo-3,6,9,12-tetraazapentadecan-15-oic acid, trifluoroacetate salt

Synonyms: Ac-Trp-Glu-Ala-Asp-pNA, Ac-Trp-Glu-Ala-Asp-p-nitroanilide, Caspase-1 and Caspase-4 Chromogenic Substrate

MF: C₃₁H₃₅N₇O₁₁ • XCF₃COOH

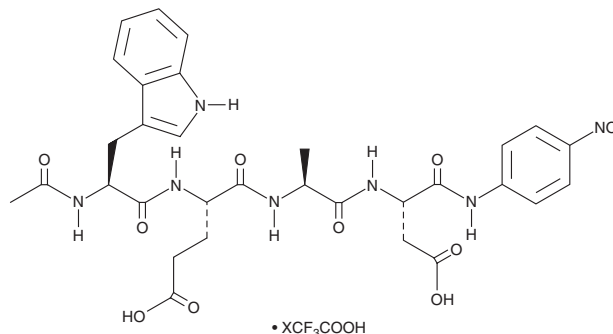
FW: 681.7

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

Description

Ac-WEAD-pNA is a colorimetric substrate for caspase-1 and caspase-4.¹ Caspase-1 and caspase-4 preferentially bind to and cleave the Trp-Glu-Ala-Asp (WEAD) peptide sequence to release p-nitroanilide (pNA), which can be quantified by colorimetric detection at 405 nm as a measure of caspase-1 and caspase-4 activity.

Reference

1. Lee, D., Adams, J.L., Brandt, M., *et al.* A substrate combinatorial array for caspases. *Bioorg. Med. Chem. Lett.* 9(12), 1667-1672 (1999).

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, 11/03/2022

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