

PRODUCT INFORMATION



Ac-WEHD-AFC (trifluoroacetate salt)

Item No. 14992

Formal Name: N-acetyl-L-tryptophyl-L- α -glutamyl-L-histidyl-N-[2-oxo-4-(trifluoromethyl)-2H-1-benzopyran-7-yl]-L- α -asparagine, trifluoroacetate salt

Synonyms: N-Acetyl-Trp-Glu-His-Asp-7-amino-4-Trifluoromethylcoumarin, Caspase-1 Substrate (Fluorogenic), Caspase-5 Substrate (Fluorogenic)

MF: C₃₈H₃₇F₃N₈O₁₁ • XCF₃COOH

FW: 838.7

Purity: \geq 95%

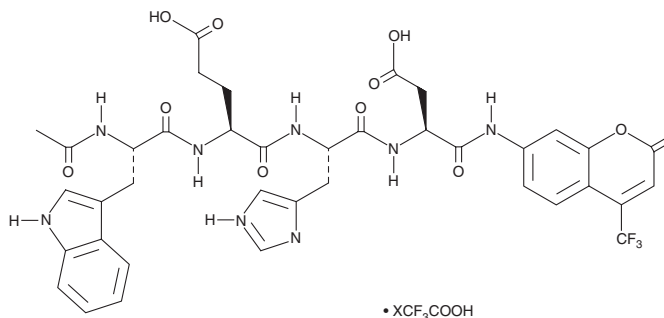
UV/Vis.: λ_{\max} : 210, 338 nm

Ex./Em. Max: 400/505 nm

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-WEHD-AFC (trifluoroacetate salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the Ac-WEHD-AFC (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Ac-WEHD-AFC (trifluoroacetate salt) is soluble in organic solvents such as DMSO. It is also soluble in water. The solubility of Ac-WEHD-AFC (trifluoroacetate salt) in DMSO and water is approximately 10 and 1 mg/ml, respectively. We do not recommend storing the aqueous solution for more than one day.

Description

Ac-WEHD-AFC is a fluorogenic substrate that can be cleaved by group I caspases (caspase-1, -4, and -5).¹ Caspase activity can be quantified by fluorescent detection of free AFC (also known as 7-amino-4-trifluoromethylcoumarin), which is excited at 400 nm and emits at 505 nm.

Reference

1. Park, K., Kuechle, M.K., Choe, Y., *et al.* Expression and characterization of constitutively active human caspase-14. *Biochem. Biophys. Res. Commun.* **347**(4), 941-948 (2006).

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, 10/18/2022

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