

# PRODUCT INFORMATION



## Ac-IETD-AFC (trifluoroacetate salt)

Item No. 17480

**Formal Name:** N-acetyl-L-isoleucyl-L- $\alpha$ -glutamyl-L-threonyl-N-[2-oxo-4-(trifluoromethyl)-2H-1-benzopyran-7-yl]-L- $\alpha$ -asparagine, trifluoroacetate salt

**Synonym:** N-Acetyl-Ile-Glu-Thr-Asp-7-amino-4-Trifluoromethylcoumarin

**MF:** C<sub>31</sub>H<sub>38</sub>F<sub>3</sub>N<sub>5</sub>O<sub>12</sub> • XCF<sub>3</sub>COOH

**FW:** 729.7

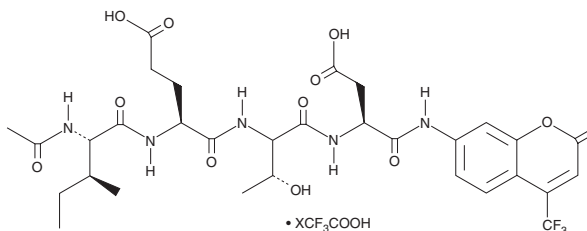
**Purity:**  $\geq 95\%$

**UV/Vis.:**  $\lambda_{\max}$ : 232, 340 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-IETD-AFC (trifluoroacetate salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the Ac-IETD-AFC (trifluoroacetate salt) in the solvent of choice. Ac-IETD-AFC (trifluoroacetate salt) is soluble in organic solvents such as DMSO, which should be purged with an inert gas, at a concentration of approximately 10 mg/ml.

Ac-IETD-AFC (trifluoroacetate salt) is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

### Description

Ac-IETD-AFC is a fluorogenic substrate that can be cleaved by caspase-8 and related enzymes that recognize the amino acid sequence IETD, including caspase-3 processing enzyme, caspase-10, and granzyme B.<sup>1</sup> 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. Reszka, P., Schulz, R., Methling, K., *et al.* Synthesis, enzymatic evaluation, and docking studies of fluorogenic caspase 8 tetrapeptide substrates. *ChemMedChem* 5, 103-117 (2010).

#### 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, 11/17/2022

#### 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