

# PRODUCT INFORMATION



## Ac-IETD-AMC (trifluoroacetate salt)

Item No. 43014

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

**Synonym:** Ac-Ile-Glu-Thr-Asp-AMC

**MF:** C<sub>31</sub>H<sub>41</sub>N<sub>5</sub>O<sub>12</sub> • XCF<sub>3</sub>COOH

**FW:** 675.7

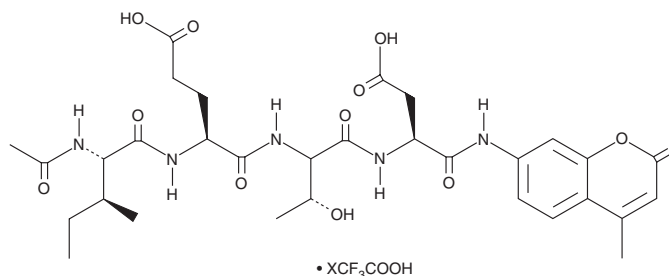
**Purity:**  $\geq$ 98%

**Supplied as:** A solid

**Ex./Em. Max:** 340-360/440-460 nm

**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-AMC (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the Ac-IETD-AMC (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Ac-IETD-AMC (trifluoroacetate salt) is soluble ( $\geq$ 10 mg/ml) in DMSO.

### Description

Ac-IETD-AMC is a fluorogenic substrate for caspase-8.<sup>1</sup> Upon enzymatic cleavage by caspase-8, 7-amino-4-methylcoumarin (AMC) is released and its fluorescence can be used to quantify caspase-8 activity. AMC displays excitation/emission maxima of 340-360/440-460 nm, respectively.

### Reference

1. Frost, V., Al-Mehairi, S., and Sinclair, A.J. Exploitation of a non-apoptotic caspase to regulate the abundance of the cdk1 p27(KIP1) in transformed lymphoid cells. *Oncogene* **20(22)**, 2737-2748 (2001).

#### 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, 05/12/2025

#### 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