

PRODUCT INFORMATION



Ac-IEPD-AFC

Item No. 42179

CAS Registry No.: 1135417-31-0
Formal Name: (S)-4-((2S,3S)-2-acetamido-3-methylpentanamido)-5-(((S)-3-carboxy-1-oxo-1-((2-oxo-4-(trifluoromethyl)-2H-chromen-7-yl)amino)propan-2-yl)carbamoyl)pyrrolidin-1-yl)-5-oxopentanoic acid

Synonym: Ac-Ile-Glu-Pro-Asp-AFC

MF: C₃₂H₃₈F₃N₅O₁₁

FW: 725.7

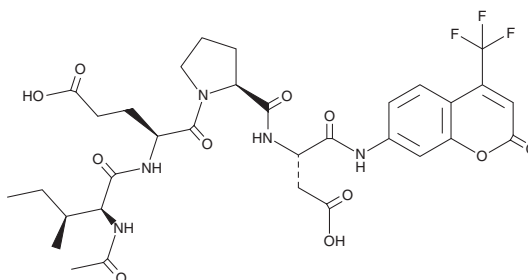
Purity: ≥98%

Ex./Em. Max: 400/505 nm

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-IEPD-AFC is supplied as a solid. A stock solution may be made by dissolving the Ac-IEPD-AFC in the solvent of choice, which should be purged with an inert gas. Ac-IEPD-AFC is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

Ac-IEPD-AFC is a fluorogenic substrate for granzyme B.¹ Upon enzymatic cleavage by granzyme B, 7-amino-4-trifluoromethylcoumarin (AFC) is released and its fluorescence can be used to quantify granzyme B activity. AFC displays excitation/emission maxima of 400/505 nm, respectively.

Reference

1. Briones, J.C., Espulgar, W.V., Koyama, S., et al. A microfluidic platform for single cell fluorometric granzyme B profiling. *Theranostics* **10**(1), 123-132 (2020).

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, 11/11/2024

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