

PRODUCT INFORMATION



Ac-LETD-AFC

Item No. 14991

CAS Registry No.: 210345-02-1

Formal Name: N-acetyl-L-leucyl-L- α -glutamyl-L-threonyl-N-[2-oxo-4-(trifluoromethyl)-2H-1-benzopyran-7-yl]-L- α -asparagine
Synonyms: N-Acetyl-Leu-Glu-Thr-Asp-7-amino-4-Trifluoromethylcoumarin, Caspase-8 Substrate (Fluorogenic)

MF: $C_{31}H_{38}F_3N_5O_{12}$

FW: 729.7

Purity: $\geq 95\%$

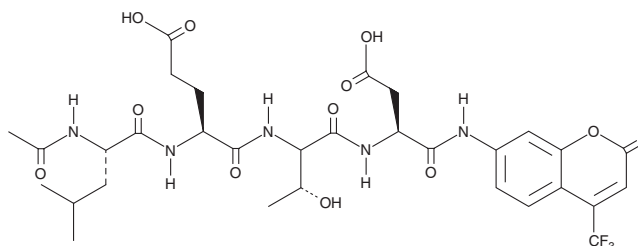
UV/Vis.: λ_{max} : 206, 339 nm

Ex./Em. Max: 400/505 nm

Supplied as: A crystalline 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-LETD-AFC is supplied as a crystalline solid. A stock solution may be made by dissolving the Ac-LETD-AFC in the solvent of choice. Ac-LETD-AFC is soluble in the organic solvent DMSO, which should be purged with an inert gas. The solubility of Ac-LETD-AFC in DMSO is approximately 10 mg/ml.

Description

Ac-LETD-AFC is a fluorogenic substrate that can be cleaved specifically by caspase-8. 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.

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.

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