

PRODUCT INFORMATION

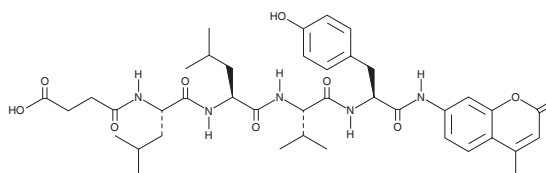


Suc-Leu-Leu-Val-Tyr-AMC

Item No. 10008119

CAS Registry No.: 94367-21-2
Formal Name: N-(3-carboxy-1-oxopropyl)-L-leucyl-L-leucyl-L-valyl-N-(4-methyl-2-oxo-2H-1-benzopyran-7-yl)-L-tyrosinamide

Synonym: Suc-LLVY-AMC
MF: C₄₀H₅₃N₅O₁₀
FW: 763.9
Purity: ≥98%
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 330 nm
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Suc-Leu-Leu-Val-Tyr-AMC is supplied as a crystalline solid. A stock solution may be made by dissolving the Suc-Leu-Leu-Val-Tyr-AMC in the solvent of choice, which should be purged with an inert gas. Suc-Leu-Leu-Val-Tyr-AMC is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of Suc-Leu-Leu-Val-Tyr-AMC in ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Suc-Leu-Leu-Val-Tyr-AMC is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, Suc-Leu-Leu-Val-Tyr-AMC should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Suc-Leu-Leu-Val-Tyr-AMC has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Suc-Leu-Leu-Val-Tyr-AMC is a fluorescent substrate for the 20S proteasome, other chymotrypsin-like proteases, and calpains (excitation max: 360 nm; emission max: 460 nm).^{1,2}

References

1. Geier, E., Pfeifer, G., Wilm, M., *et al.* A giant protease with potential to substitute for some functions of the proteasome. *Science* **283**(5404), 978-981 (1999).
2. Sasaki, T., Kikuchi, T., Yumoto, N., *et al.* Comparative specificity and kinetic studies on porcine calpain I and calpain II with naturally occurring peptides and synthetic fluorogenic substrates. *J. Biol. Chem.* **259**(20), 12489-12494 (1984).

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, 11/30/2022

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