Suc-Leu-Leu-Val-Tyr-AMC

Proteasome substrate

Fluorogenic substrate for 20S proteasome (Prod. No. BML-PW8720) and other chymotrypsin-like proteases, as well as calpains. Ex.: 380nm Em.: 460nm.

Citations: 105

View Online »

Ordering Information

Order Online »

BML-P802-0005

5mg

Manuals, SDS & CofA

View Online »

- Fluorogenic substrate of calpains and 20S proteasome
- Used to study the activity of calpains and proteosomal degradation

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Handling Protect from light.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Suc-LLVY-AMC, Proteasome substrate (fluorogenic),

Calpain substrate (fluorogenic)

CAS 94367-21-2

Formula $C_{40} H_{53} N_5 O_{10}$

Formulation Lyophilized.

Identity Determined by ESI-MS.

MW 763.9

Purity ≥97% (HPLC)

Sequence N-Succinyl-Leu-Leu-Val-Tyr-AMC (AMC=7-amino-4-

methylcoumarin)

Solubility Soluble in DMSO (at least 5mM).

Last modified: May 29, 2024

