Z-Leu-Leu-Leu-vinyl sulfone

Inhibits the trypsin-like, chymotrypsin-like and caspase-like (PGPH) activity in isolated proteasomes and living cells by covalent modification of the amino-terminal threonine of the catalytically active β subunits. The vinyl sulfone acts as a Michael acceptor for soft nucleophiles such as thiols, leading to the formation of a covalent bond, while being resistant to attack by free thiols, which can cause inactivation of other classes of protease inhibitors.

Ordering Information

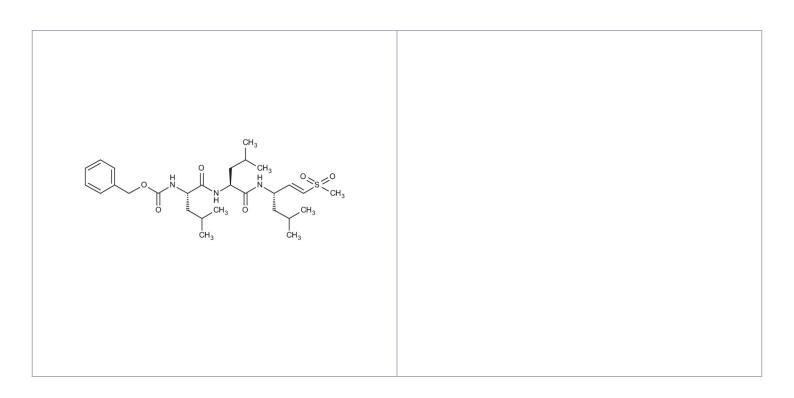
Order Online »

BML-ZW9170-0500

500µg

Manuals, SDS & CofA

View Online »



Handling & Storage

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

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

MW ~551.3 Da

Purity ≥99%

Sequence Z-(Leu)₃-vinyl sulfone

Last modified: May 29, 2024

