

Ada-(Ahx)₃-(Leu)₃-vinyl sulfone

A new class of proteasome inhibitors, considerably extended in comparison with the commonly used fluorescent substrates and peptide-based inhibitors, has been prepared. Evaluation of these new compounds revealed a set of highly potent proteasome inhibitors that target all individual active subunits with comparable affinity, unlike the other inhibitors described to date. These compounds are more potent in living cells than their shorter peptide vinyl sulphone counterparts. Modification of the most active compound, adamantaneacetyl-(6-aminohexanoyl)3-(leucyl)3-vinylmethyl-sulphone itself capable of proteasome inhibition in living cells, afforded a new set of radio- and affinity labels.

Citations: 1

[View Online »](#)

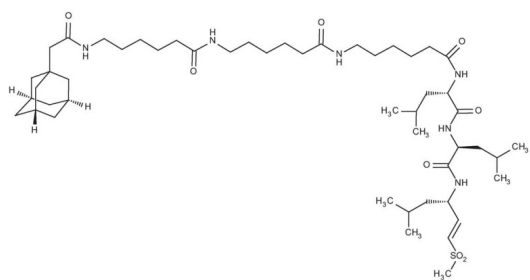
Ordering Information

[Order Online »](#)

BML-AW9155-0100	100µg
------------------------	-------

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

Appearance White lyophilized powder.

Formula $C_{50}H_{88}N_6O_8S$

MW 933.1

Purity ≥94% (HPLC)

Sequence Ada-(Ahx)₃-(Leu)₃-vinyl sulfone

Solubility Soluble in formic acid (1 mg/ml).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)