

Proteasome inhibitor I (aldehyde)

Cell permeable inhibitor of the 26S proteasome (MCP; multicatalytic proteinase complex) in HT4 cells. Causes accumulation of ubiquitinated proteins in neuronal cells. Prevents the activation of NF- κ B in response to TNF- α or okadaic acid (Prod. No. ALX-350-003) through inhibition of I κ B α degradation.

Citations: 4

[View Online »](#)

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 Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Z-IE(OtBu)AL-CHO, Z-Ile-Glu(OtBu)-Ala-Leu-CHO, PSI

Appearance White powder.

Formula $C_{32}H_{50}N_4O_8$

MW 618.8

Purity ≥95% (TLC)

Sequence Z-Ile-Glu(OtBu)-Ala-Leu-CHO

Solubility Soluble in 100% ethanol (30mg/ml), DMSO (30mg/ml) or methanol (50mg/ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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