Ac-Ala-Pro-Nle-Asp-OH

Highly specific inhibitor of the caspase-like peptidic activity of the 20S proteasome.

Citations: 1

View Online »

Ordering Information

Order Online »

BML-AW9485-0100

100µg

Manuals, SDS & CofA

View Online »

Handling & Storage

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

use solutions immediately and do not store.

Handling Do not add buffers containing free amino groups as these will react with the aldehyde

moiety thus inactivating the compound.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance Glassy lyophilized solid.

Formula $C_{20}H_{32}N_4O_8$

MW 456.5

Purity ≥95% (Reverse-Phase HPLC)

Sequence Ac-Ala-Pro-Nle-Asp-OH

Solubility Soluble in water:acetonitrile containing 0.1% TFA (3:1) at a

concentration of 1mg/mL.

Last modified: May 29, 2024