# **Z-DEVD-pNA**

### **Caspase substrate**

Chromogenic substrate for caspase-3 with increased cell permeability. Also substrate for caspase-1, -4 and -7.  $\lambda_{max}$  of pNA is 400nm. Similar to Ac-DEVD-pNA except the N-terminus is blocked by a benzyloxycarbonyl (Z) moiety which may increase cell permeability.

Citations: 5

View Online »

**Ordering Information** 

Order Online »

ALX-260-048-M001

1mg

Manuals, SDS & CofA

**View Online** »

# **Handling & Storage**

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

**Handling** Keep cool and dry.

Long Term Storage -20°C

**Shipping** Ambient Temperature

## Regulatory Status RUO - Research Use Only

#### **Product Details**

Alternative Name Caspase-3 substrate (chromogenic)

**Appearance** White powder.

Formula  $C_{32}H_{38}N_6O_{14}$ 

Formulation Lyophilized.

**MW** 730.7

Peptide Content 70-90%

Purity ≥99% (HPLC)

**Sequence** Z-Asp-Glu-Val-Asp-pNa (pNA = p-Nitroaniline)

**Soluble** in DMSO or acetonitrile.

Last modified: May 29, 2024