Ac-VDVAD-AFC

Caspase-2 substrate

Fluorogenic substrate for caspase-2; similar to Ac-VDVAD-AMC but the AFC fluorophore has a greater Stokes' shift upon cleavage. VDVAD has been found to be a preferred cleavage site for caspase-2 (Ich-1L). Increases in Ac-VDVAD-AFC cleavage correlated with losses in procaspase-2 (conversion to active caspase-2) in human neuroblastoma lines induced to apoptosis with C2-ceramide and NO.

Ordering Information

Order Online »

ALX-260-112-M005	5mg
ALX-260-112-M010	10mg

Manuals, SDS & CofA

View Online »

Handling & Storage

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

Handling Warm up to room temperature before opening. Keep cool and dry.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Caspase-2 substrate (fluorogenic)

Appearance White to off-white powder.

Formula $C_{33}H_{41}F_3N_6O_{12}$

MW 770.7

Peptide Content 75-95%

Purity ≥98% (HPLC)

Sequence Ac-Val-Asp-Val-Ala-Asp-AFC (AFC=7-Amino-4-

trifluoromethylcoumarin)

Solubility Soluble in DMSO.

Technical Info / Product NotesAFC has an excitation maximum of 400nm and an

emission maximum of 505nm.

Last modified: May 29, 2024

