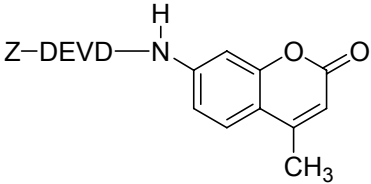


# Caspase-3 Substrate Ac-DEVD-AMC

7-Amino-4-methylcoumarin, N-CBZ-L-aspartyl-L-glutamyl-L-valyl-L-aspartic acid amide (Z-DEVD-MCA)

## Instruction Manual

Catalog Number	PK-CA707-10202
Description	Z-DEVD-AMC is a fluorogenic substrate for caspase 3 and can be used to continuously measure the activity of caspase-3 in cell extracts, using a fluorometer or fluorescence microplate reader.
Quantity	5 mg
Excitation / Emission Maxima	$\lambda_{exc}/\lambda_{em} = 330/390$ nm (of substrate, weak fluorescence); $\lambda_{exc}/\lambda_{em} = 342/441$ nm (of end product AMC, strong fluorescence)
Molecular Structure	
Molecular Weight / Molecular Formula	675.74 Da; C <sub>30</sub> H <sub>37</sub> N <sub>5</sub> O <sub>13</sub>
Purity	>95% (as determined by TLC)
Appearance / Formulation / Solubility	Off-white solid; soluble in DMSO.
Storage & Stability	Store desiccated at -20°C. Protect from light.
Applications	For use in caspase-3 activity assays.
References	1) Anal Biochem 292, 313 (2001) 2) Biochemistry 39, 16056(2000) 3) Br J Pharmacol 129, 1673(2000) 4) J Biol Chem 275, 21094(2000)
Caution	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.