



Product Data Sheet

Product Name:	Caspase 3 (Apopain) Substrate 1f, fluorogenic	
Catalog Number:	AS-25273-5 (5 mg)	Lot Number: See label on vial
Sequence:	Ac-DEVD-AFC (1-letter code)	
Molecular Weight:	729.5	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	

Peptide Reconstitution: Caspase 3 peptide is freely soluble in basic buffer.

Storage: Caspase 3 peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at -20°C or lower. Reconstituted peptide can be aliquoted and stored at -20°C or lower.

Description: AFC (7-amino-4-trifluoromethylcoumarin)-derived caspase substrates are widely used for the fluorimetric detection of various caspase activities. Cleavage of AFC peptides by caspases generate strongly fluorescent AFC that is monitored fluorimetrically at 500-510 nm with excitation of 370-390 nm. Caspase-3 substrate with $k_{\text{cat}}/K_{\text{m}} = 1.3 \times 10^6 \text{ M}^{-1}\text{s}^{-1}$, Caspase-7 substrate with $k_{\text{cat}}/K_{\text{m}} = 1.3 \times 10^4 \text{ M}^{-1}\text{s}^{-1}$. Ref: Gurtu V, et al. *Anal Biochem* **251**, 98 (1997); Martins LM, et al. *J Biol Chem* **272**, 7421 (1997); Deveraux QL, et al. *Nature* **388**, 300 (1997); Xiang J, et al. *Proc Natl Acad Sci USA* **93**, 14559 (1997).

Related Products:

Name	Cat #	Size
SensoLyte® Homogeneous AFC Caspase-3/7 Assay Kit 500 assays (96-well plate)	AS-71114	1 kit
Caspase 3 (Apopain) Substrate 1, chromogenic	AS-25263-5	5 mg
Caspase 3 (Apopain) Substrate 1m, fluorogenic	AS-25262-5	5 mg

For Research Use Only