

Ac-DEVD-AMC

Ac-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. For real-time detection of caspase-3 activity in intact cells, see our novel [NucView@488 Caspase-3 Assay Kit](#).



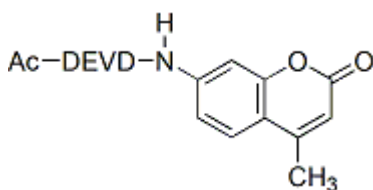
Product attributes

CAS number	169332-61-0
Apoptosis/viability marker	Caspase
For live or fixed cells	Cell lysis required
Detection method/readout	Microplate reader (fluorescence)
Assay type/options	Endpoint assay
Substrate specificity	Caspases
Colors	Blue
Excitation/Emission	342/441 nm (end product)

Product Description

Ac-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. For real-time detection of caspase-3 activity in intact cells, see our novel [NucView@488 Caspase-3 Assay Kit](#).

- Ex/Em of substrate = 330/390 nm (weak fluorescence)
- Ex/Em of end product (AMC) = 342/441 nm (strong fluorescence)
- Off-white solid soluble in DMSO
- C₃₀H₃₇N₅O₁₃
- MW: 675.74



Biotium Catalog #10202

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).

This datasheet was generated on June 19, 2026 at 08:25:29 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/ac-devd-amc-7-amino-4-methylcoumarin-n-acetyl-l-aspartyl-l-glutamyl-l-valyl-l-aspartic-acid-amide/>