

Ac-DEVD-R110

A fluorogenic substrate for caspase-3 that can be used to continuously measure the activity of caspase-3 in cell extracts, using a fluorometer or fluorescence microplate reader.

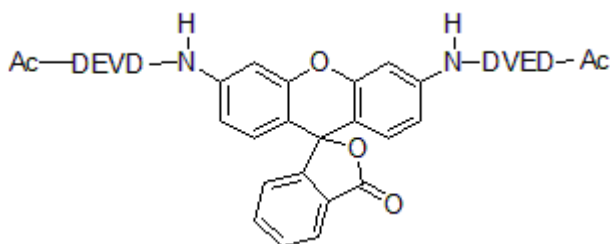


Product Description

Ac-DEVD-R110 (Rhodamine 110, bis-(N-Ac-L-aspartyl-L-glutamyl-L-valyl-1-aspartic acid amide)) 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. Ac-DEVD-R110 contains two DEVD tetrapeptides and is hydrolyzed by caspase-3/7 in two successive steps to release the green fluorescent dye rhodamine 110 (R110).

For caspase-3 kits using Ac-DEVD-R110, please see our Caspase-3 DEVD-R110 Fluorometric and Colorimetric Assay Kit ([30008](#)), and our Caspase-3 DEVD-R110 Fluorometric HTS Assay Kit ([30009](#)). For real-time detection of caspase-3 activity in intact cells, see our novel [NucView@488 Caspase-3 Assay Kit](#).

- Abs/Em of end product (R110) = 496/520 nm (strong fluorescence)
- Off-white solid soluble in DMSO
- C₆₀H₇₀N₁₀O₂₅
- MW: 1331



Product attributes

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	Green
Excitation/Emission	496/520 nm

References

1. Cell Death Differ 6(2), 99(1999) [doi:10.1038/sj.cdd.4400476](https://doi.org/10.1038/sj.cdd.4400476)
2. J Biol Chem 275, 288(2000) [doi:10.1074/jbc.M112.447540](https://doi.org/10.1074/jbc.M112.447540)
3. Biochemistry 38(42), 13906(1999) [doi:10.1021/bi9913395](https://doi.org/10.1021/bi9913395)

This datasheet was generated on June 15, 2026 at 08:21:04 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/ac-devd-r110/>