

Z-AAD-R110

Z-AAD-R110 is a substrate for granzyme B. Granzyme B from T lymphocytes and natural killer cells induces apoptosis in target cells by activating caspases and causing mitochondrial cytochrome C release.



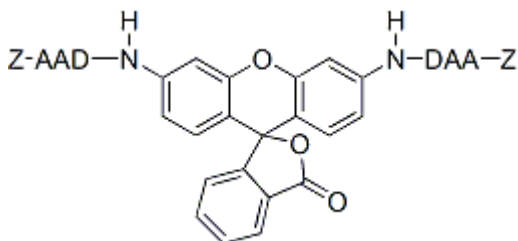
Product attributes

Excitation/Emission	496/520 nm (end product)
---------------------	--------------------------

Product Description

Z-AAD-R110 is a substrate for granzyme B. Granzyme B from T lymphocytes and natural killer cells induces apoptosis in target cells by activating caspases and causing mitochondrial cytochrome C release.

- Full name: Rhodamine 110, bis-(N-CBZ-L-alanyl-L-alanyl-L-aspartic acid amide)
- $\lambda_{Ex}/\lambda_{Em}$ of end product (R110) = 496/520 nm
- Off-white to pink solid soluble in DMSO
- Store at -20°C and protect from light
- $\text{C}_{56}\text{H}_{56}\text{N}_8\text{O}_{17}$
- MW: 1113.1



References

1. J Biol Chem 276, 6974 (2001).

This datasheet was generated on June 10, 2026 at 03:43:43 AM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/z-aad-r110-rhodamine-110-bis-n-cbz-l-alanyl-l-alanyl-l-aspartic-acid-amide/>