

Ac-IETD-R110

Z-IETD-R110 is a substrate for caspase-8, which is involved at the early stage of apoptosis, acting as an initiator of the caspase activation cascade.



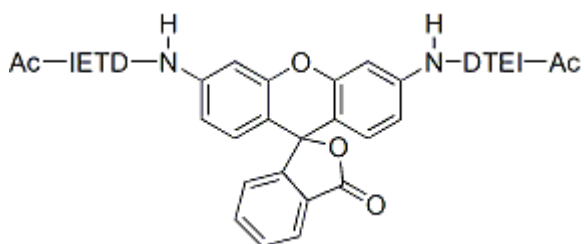
Product attributes

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

Product Description

Z-IETD-R110 is a substrate for caspase-8, which is involved at the early stage of apoptosis, acting as an initiator of the caspase activation cascade.

- Full name: Rhodamine 110, bis-(N-acetyl-L-isoleucyl-L-glutamyl-L-threonyl-L-aspartic acid amide
- $\lambda_{Ex}\lambda_{Em}$ of end product (R110) = 496/520 nm
- Off-white to pink solid soluble in DMSO
- $C_{68}H_{82}N_{10}O_{24}$
- MW: 1422



This datasheet was generated on June 10, 2026 at 02:20:36 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/ac-ietd-r110-rhodamine-110-bis-n-acetyl-l-isoleucyl-l-glutamyl-l-threonyl-l-aspartic-acid-amide/>