



Glowing products for science

Caspase-8 IETD-R110 Fluorometric HTS Assay Kit

This item is discontinued.

Product Description

This item is discontinued.



Call us : [800-304-5357](tel:800-304-5357)

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, High-throughput assay, Homogeneous assay
Substrate specificity	Caspases
Colors	Green
Excitation/Emission	496/520 nm (end product)
Storage Conditions	Store at -10 to -35 °C

This datasheet was generated on January 7, 2026 at 10:23:38 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/caspase-8-ietd-r110-fluorometric-hts-assay-kit/>