



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

<b>Apoptosis/viability marker</b>	Caspase
<b>For live or fixed cells</b>	Cell lysis required
<b>Detection method/readout</b>	Microplate reader (fluorescence)
<b>Assay type/options</b>	Endpoint assay, High-throughput assay, Homogeneous assay
<b>Substrate specificity</b>	Caspases
<b>Colors</b>	Green
<b>Excitation/Emission</b>	496/520 nm (end product)
<b>Storage Conditions</b>	Store at -10 to -35 °C

This datasheet was generated on June 8, 2026 at 07:48:36 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/>