



Glowing products for science

Caspase-8 IETD-R110 Fluorometric and Colorimetric Assay Kit

This item is discontinued.

Product Description

This item is discontinued.

References

1. Neurochem Res (2009) Aug;34(8):1410-6. doi: [10.1007/s11064-009-9927-6](https://doi.org/10.1007/s11064-009-9927-6)
2. J Psychiatr Res (2009) May;43(8):754-62 doi: [10.1016/j.jpsychires.2008.10.002](https://doi.org/10.1016/j.jpsychires.2008.10.002)
3. Eur J Pharmacol (2009) May 21;610(1-3):99-105 doi: [10.1016/j.ejphar.2009.03.046](https://doi.org/10.1016/j.ejphar.2009.03.046)
4. Fundam Clin Pharmacol (2010) Aug;24(4):445-50 doi: [10.1111/j.1472-8206.2009.00789.x](https://doi.org/10.1111/j.1472-8206.2009.00789.x)
5. Asian Pac J Cancer Prev (2018) 19(4): 917–922 doi: [10.22034/APJCP.2018.19.4.917](https://doi.org/10.22034/APJCP.2018.19.4.917)

This datasheet was generated on August 15, 2026 at 09:47:38 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/caspase-8-ietd-r110-fluorometric-and-colorimetric-assay-kit/>



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 (absorbance), Microplate reader (fluorescence)
Assay type/options	Endpoint assay
Substrate specificity	Caspases
Colors	Green
Excitation/Emission	496/520 nm (end product)
Storage Conditions	Store at -10 to -35 °C