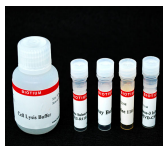


## Caspase-3 DEVD-R110 Fluorometric and Colorimetric Assay Kit

This item is discontinued.



### 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 (absorbance), Microplate reader (fluorescence)
<b>Assay type/options</b>	Endpoint 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

### Product Description

This item is discontinued.

### References

1. FASEB J (2008) Jul;22(7):2243-52. doi: [10.1096/fj.07-099234](https://doi.org/10.1096/fj.07-099234)
2. Plant Biol (2008) Nov;10(6):732-45 doi: [10.1111/j.1438-8677.2008.00078.x](https://doi.org/10.1111/j.1438-8677.2008.00078.x).
3. BMC Immunol (2009) 10, 29 doi:[10.1186/1471-2172-10-29](https://doi.org/10.1186/1471-2172-10-29)
4. J Plant Physiol (2012) Oct 15;169(15):1489-500 doi: [10.1016/j.jplph.2012.06.002](https://doi.org/10.1016/j.jplph.2012.06.002)
5. Growth Factors (2012) Dec;30(6):385-93 doi: [10.3109/08977194.2012.734506](https://doi.org/10.3109/08977194.2012.734506)
6. Clin Exp Immunol (2012) Aug;169(2):156-63 doi: [10.1111/j.1365-2249.2012.04607.x](https://doi.org/10.1111/j.1365-2249.2012.04607.x).
7. J Cell Physiol (2013) Jul;228(7):1568-76 doi: [10.1002/jcp.24318](https://doi.org/10.1002/jcp.24318).
8. Eur J Pharmacol (2013) Jan 5;698(1-3):87-94 doi: [10.1016/j.ejphar.2012.10.007](https://doi.org/10.1016/j.ejphar.2012.10.007).
9. Eur J Pharmacol (2014) Oct 5;740:683-9 doi: [10.1016/j.ejphar.2014.06.009](https://doi.org/10.1016/j.ejphar.2014.06.009).
10. Nutr Cancer (2020) Apr 9:1-11 doi: [10.1080/01635581.2020.1751217](https://doi.org/10.1080/01635581.2020.1751217)

This datasheet was generated on June 15, 2026 at 09:49:34 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/caspase-3-devd-r110-fluorometric-colorimetric-assay-kit/>