Dual Apoptosis Assay with NucView® 488 Caspase-3 Substrate & Annexin V

These kits provides two apoptosis markers, our novel NucView® 488 Caspase-3 Substrate and Annexin V conjugate for detecting caspase-3 activation and phosphatidylserine (PS) translocation.

Product Description

These kits provides two apoptosis markers, our novel NucView® 488 Caspase-3 Substrate and fluorescent CF® dye Annexin V for detecting two important apoptosis events, caspase-3 activation and phosphatidylserine (PS) translocation.

The rate of apoptosis varies from cell to cell even within the same population. As a result, various apoptotic events or markers accompanying the apoptotic process also occur differently among cells. Thus, it is important to be able to detect these apoptotic events on an individual cell basis. Traditionally, caspase activity has been detected either using a membrane-impermeable fluorogenic enzyme substrate such as DEVD-R110, or a fluorescently-labeled inhibitor such as a FLICA reagent. In the former case, cell lysis is required, thus precluding the detection of caspase activity in live cells. In addition, such caspase assays measure only the average caspase activity of a highly heterogeneous cell population at a given time. In the latter case, although a FLICA reagent can enter live cells to detect caspase activity, only the initial fluorescent signal following the application of the reagent can truly reflect the enzyme activity or the state of the apoptotic cells because any detected signal after the initial "snapshot" will need to consider the potential interference of the inhibitor to the enzyme and the apoptotic cell itself.

Unlike conventional caspase assays, NucView® 488 Caspase-3 substrate detects caspase-3 activity within individual whole cells in a non-interfering manner. The substrate consists of a fluorogenic DNA dye and a DEVD substrate moiety specific for caspase-3. The substrate, which is both non-fluorescent and nonfunctional as a DNA dye, rapidly crosses cell membranes to enter the cytoplasm, where it is cleaved by caspase-3 to form a high-affinity DNA dye that stains the nucleus bright green. Thus, the NucView® 488 caspase-3 substrate is bi-functional, allowing detection of caspase-3 activity and visualization of apoptotic nuclear morphology

The kit contains reagents sufficient for 50 flow cytometry assays (200 uL assay volume). The number of fluorescence microscopy assays that can be performed with the kit may vary based on the size of culture vessel and staining volume used.

To learn about the advantages of monitoring apoptosis using NucView® caspase-3 substrates, visit the NucView® Technology Page.

NucView is a registered trademark of Biotium. NucView enzyme substrate technology is covered by US patents. Texas Red is a registered trademark of Thermo Fisher Scientific.

References

- 1. Reproductive BioMedicine Online (2008) 16(5), 657-663. www.rbmonline.com/Article/3185
- 2. Curr Trends Biotech Pharm (2009) 3(3), 278-286. ISSN 0973-8916
- 3. Malaysian Journal of Chemistry (2009) Vol 11(1), 136-142.
- 4. Malaysian Journal of Chemistry (2009) 11(1), 143 148.
- 5. Pharmacognosy Res (2010) 2(2), 113-119. doi: 10.4103/0974-8490.62949
- 6. Mol Cancer Res (2013) 11(4), 405-17. DOI: 10.1158/1541-7786.MCR-12-0551
- Stem Cells (2013) 31, 1121-1135. doi: 10.1002/stem.1368
- 8. Stem Cells (2013) 31, 2374-2387. https://doi.org/10.1002/stem.1509
- 9. Journal of Cell Science (2014) 127, 1738-1750. doi: 10.1242/jcs.138214
- 10. PLoS Genet (2014) 10(10), e1004594. doi:10.1371/journal.pgen.1004594
- 11. Cancer Gene Therapy (2016) Cancer Gene Ther 23(6), 178-87. doi:10.1038/cgt.2016.18
- 12. Current Cancer Drug Targets (2018) 18, 1-9. DOI:10.2174/1568009617666171114144236 13. Sci Rep (2018) 8, 13192. DOI:10.1038/s41598-018-31575-3

Find a list of NucView® references and a list of validated cell lines under Supporting Documents.

Apoptosis/viability marker For live or fixed cells For live/intact cells

Product attributes

Call us: 800-304-5357

Fluorescence microscopy, Live cell imaging, Flow cytome Detection method/readout Assay type/options Endpoint assay

Phosphatidylserine/Annexin V, Caspase

Email: btinfo@biotium.com

Fixation options Fix after staining (formaldehyde)

Green/Red, Green/Far-red Product origin Annexin V (human); recombinant, produced in E. coli

Storage Conditions Store at 2 to 8 °C, Do not freeze, Protect from light

This datasheet was generated on November 18, 2025 at 03:08:52 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/dual-apoptosis-assay-with-nucview-488-caspase-3-substrate-and-annexin-v/