## **Resazurin Cell Viability Assay Kit** (alamarBlue™)

A sensitive, safe and cost-effective high-throughput homogeneous assay of cell viability by fluorescence or absorbance measurement.

## **Product Description**

Resazurin Cell Viability Assay offers a simple, rapid, reliable, sensitive, safe and cost-effective measurement of cell viability. This assay has excellent performance compared to other resazurin-based cell proliferation kits such as alamarBlue®, PrestoBlue®, or CellTiter-Blue®.

- Single-step, homogeneous, high-throughput cell quantitation
- For fluorescence- or absorbance-based measurement
- Sensitive, requires as few as 80 cells for reproducible results
- Widely used in cultured mammalian cells, bacteria, protozoa, yeast, & fungi
- · Cell lysis not required, allows long-term kinetic monitoring
- Non-toxic, non-radioactive, non-carcinogenic, simple disposal
- Ready-to-use solution

Resazurin detects cell metabolic by converting from a nonfluorescent dye to the highly red fluorescent dye resorufin in response to chemical reduction of growth medium resulting from cell growth. Continued cell growth maintains a reduced environment while inhibition of growth maintains an oxidized environment. Reduction related to growth causes the redox indicator to change from the oxidized form (nonfluorescent, blue) form to the reduced form (fluorescent, red).

Reduced resazurin can be detected using fluorescence (Ex/Em 530-560/590 nm) or absorbance at 570 nm. The fluorescent or colorimetric signal is proportional to the number of living cells in the sample. Results can obtained between 1-24 hours after adding substrate. Resazurin solution is supplied ready-to-use in 25 mL (2500 assay) or 100 mL (10,000 assay) sizes.

See our full selection of Cell Viability & Apoptosis Assays.

alamarBlue is a registered trademark of Trek Biosystems. CellTiter-Blue is a registered trademark of Promega Corporation. PrestoBlue is a registered trademark of Thermo Fisher Scientific.

## References

- 1. Arterioscler, Thromb, Vasc Biol (2010) 30(10), 1940-1948. DOI:10.1161/ATVBAHA.110.205997
- 2. J Clin Endocrinol Metab, October 2010, 95(10):E181–E191 doi: 10.1210/jc.2010-0581
- 3. Nucleic Acids Res (2010) 38(22), 8131-8140. doi:10.1093/nar/gkq697
- 5. Noticle Action has (2010) 30(22), 6131–6140. <u>ubi-10.1033/hal/pkq097</u>.
  4. Scand J Clin Lab Invest. (2010) 70(7), 512-8. <u>doi: 10.3109/00365513.2010.521255</u>
  5. Tissue Eng A (2011) 17(1-2), 37-44. <u>http://doi.org/10.1089/ten.tea.2010.0188</u>
  6. Biomaterials (2011) 32(1), 128-36. <u>doi: 10.1016/j.biomaterials.2010.09.006</u>.
- 7. Food Chem (2012) 135(3), 1533-1538. https://doi.org/10.1016/j.foodchem.2012.06.058
  8. Drug Des Dev Therapy (2013) 7, 529-544. http://dx.doi.org/10.2147/DDDT.S45162
- 9. Langmuir (2013) 29(46), 14254-64. doi: 10.1021/la403533r
- 10. PLoS ONE (2013) 8(10), e76264. doi:10.1371/journal.pone.0076264
- 11. ACS Nano (2015) 9(2), 1219-1235. https://doi.org/10.1021/nn504890z
- 12. Antimicrob Agents Chemother (2016) 60(7), 4183-4196. doi:10.1128/AAC.03021-15
- 13. Curr Prot Toxicol (2016) 68, 2.24.1-2.24.15. doi: 10.1002/cptx.1
- 14. J Biol Chem (2016) 291, 21869-21879. doi: 10.1074/jbc.M115.712166
- 15. Curr Prot Stem Cell Biol (2018) 26(1), 5C.3.1-5C.3.19. DOI:10.1002/9780470151808.sc05c03s26
- 16. Int J Biol Macromol (2018) 113(1), 132-141. https://doi.org/10.1016/j.ijbiomac.2018.02.069

This datasheet was generated on November 2, 2025 at 01:40:33 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/resazurin-cell-viability-assay-kit/

## Product attributes

| Apoptosis/viability marker | Metabolic activity                                                          |
|----------------------------|-----------------------------------------------------------------------------|
| For live or fixed cells    | For live/intact cells                                                       |
| Detection method/readout   | Microplate reader (absorbance), Microplate reader (fluorescence)            |
| Assay type/options         | Endpoint assay, High-throughput assay, Homogeneous assay, Real-time imaging |
| Colors                     | Red                                                                         |
| Storage Conditions         | Store at 2 to 8 °C                                                          |
| Excitation/Emission        | 530-560/590 nm, 570 nm (for absorbance-based assay)                         |

Call us: 800-304-5357 Email: btinfo@biotium.com