

Product datasheet

Cell Viability Assay Kit (Bioluminescent) ab65314

2 References

Overview

Product name	Cell Viability Assay Kit (Bioluminescent)
Detection method	Luminescent
Sample type	Adherent cells, Suspension cells
Assay type	Direct
Sensitivity	> 10 cells/well
Range	10 cells/well - 100 cells/well
Assay time	0h 30m
Product overview	<p>Abcam's Cell Viability Assay Kit (Bioluminescent) utilizes bioluminescent detection of the ATP levels for a rapid screening of apoptosis and cell proliferation simultaneously in mammalian cells. The assay utilizes luciferase to catalyze the formation of light from ATP and luciferin, and the light can be measured using a luminometer or Beta Counter. The assay can be fully automatic for high throughput (10 seconds/sample). The high sensitivity of this assay has led to many other applications for detecting ATP production in various enzymatic reactions, as well as for detecting low level bacterial contamination in samples such as blood, milk, urine, soil, and sludge. Visit our FAQs page for tips and troubleshooting.</p>
Notes	<p>Cell death (especially apoptosis) is an energy-dependent process that requires ATP. As ATP levels fall to a point where the cell can no longer perform basic metabolic functions, the cell will die. A typical apoptotic cell exhibits a significant decrease in ATP level. Therefore, loss of ATP level in cell has been used as an indicator of cell death. In contrast, cell proliferation has been recognized by increased levels of ATP.</p>
Platform	Microplate reader

Properties

Storage instructions Store at +4°C. Please refer to protocols.

Components	200 tests
ATP (MW 551)	1 x 1mg
ATP Monitoring Enzyme (Lyophilised)	1 vial
Enzyme Reconstitution Buffer	1 x 2ml

Components	200 tests
Nucleotide Releasing Buffer	1 x 20ml

Relevance

Cell viability is a determination of living or dead cells, based on a total cell population. Cell viability assess healthy cells in a sample, with no distinction between dividing or quiescent cells. An increase in cell viability indicates cell growth, while a decrease in viability can generally be interpreted as the result of either toxic effects of compounds/agents or suboptimal culture conditions.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors