

# XTT Assay Kit ab232856

[8 References](#) [1 Image](#)

### Overview

<b>Product name</b>	XTT Assay Kit
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Adherent cells, Suspension cells
<b>Assay type</b>	Cell-based (quantitative)
<b>Product overview</b>	XTT Assay Kit ab232856 provides an easy to use tool for use with cell cultures to study changes in the number of cells and their metabolic activity. This provides insights into cell viability, cell proliferation, and cytotoxicity.

The XTT assay is based on the extracellular reduction of XTT by NADH produced in the mitochondria via trans-plasma membrane electron transport and an electron mediator.

Reduction of XTT during the assay produces a water-soluble orange-colored formazan product which dissolves directly into the culture medium, eliminating the need for an additional solubilization step (as is required in the [MTT assay](#), but not in the [MTS assay](#)).

The formazan product can be detected using an absorbance-based microplate reader.

As the reagents used in the XTT and their products are non-toxic and the assay does not involve cell lysis, the XTT assay can be used to measure multiple timepoints or for continual measurement.

XTT (2,3-bis-(2-methoxy-4-nitro-5-sulfophenyl)-2H-tetrazolium-5-carboxanilide) is a tetrazolium-based compound like the reagents used in the [MTT assay](#) and the [MTS assay](#).

Review our [cell health assay guide](#) to learn about our kits to perform a [cell viability assay](#), [cytotoxicity assay](#) or [cell proliferation assay](#).

<b>Platform</b>	Microplate reader
-----------------	-------------------

### Properties

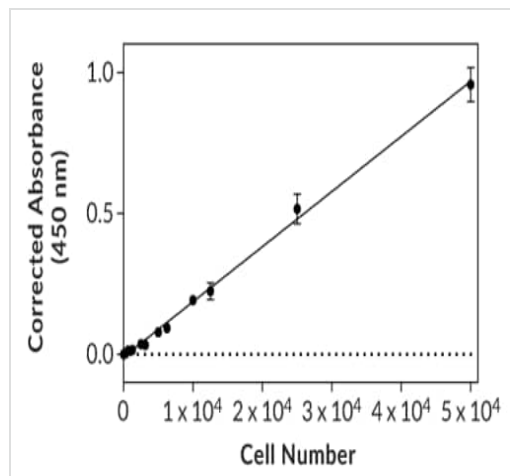
<b>Storage instructions</b>	Store at -20°C. Please refer to protocols.
-----------------------------	--

Components	1 x 96 tests
Electron Mediator Solution	1 x 600µl
XTT Developer Reagent	1 x 600µl

## Relevance

Cell proliferation is the multiplication or reproduction of cells, as a result of cell growth and cell division, resulting in the expansion of a cell population.

## Images



A typical cell titration experiment using HL-60 target cells and XTT assay kit ab232856.

A typical cell titration experiment using HL-60 target cells and XTT assay kit ab232856

**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