

SRB Assay / Sulforhodamine B Assay Kit ab235935

8 References 1 Image

Overview

Product name	SRB Assay / Sulforhodamine B Assay Kit
Detection method	Colorimetric
Sample type	Adherent cells
Assay type	Cell-based (quantitative)
Product overview	<p>SRB Assay / Sulforhodamine B Assay Kit ab235935 relies on the stoichiometric binding of SRB (Sulforhodamine B) dye to proteins under mild acidic conditions and its subsequent extraction under basic conditions. The amount of dye extracted is a proxy for cell mass and thus the number of cells in a sample.</p> <p>The measurement of the number of cells in the assay allows its use for analysis of cell viability, cytotoxicity and cell proliferation.</p> <p>In the assay, cultured cells are fixed on plates, stained with Sulforhodamine B, washed and dried, then the bound dye is solubilized and the absorbance of the dye in solution is measured at OD 565 nm.</p> <p>Review our <a href="#">cell health assay guide</a> to learn about our kits to perform a <a href="#">cell viability assay</a>, <a href="#">cytotoxicity assay</a> or <a href="#">cell proliferation assay</a>.</p>
Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K943 Sulforhodamine B Cell Cytotoxicity Assay Kit (Colorimetric). K943-1000 is the same size as the 1000 test size of ab235935.</p>
Platform	Microplate reader

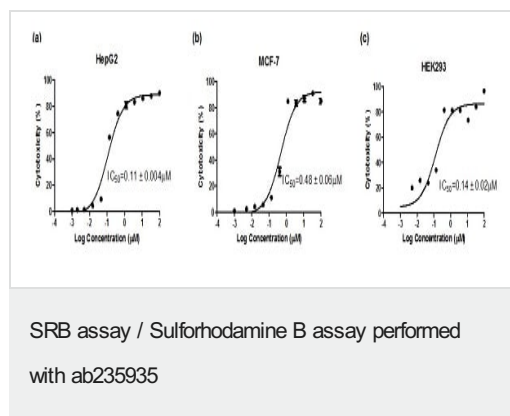
Properties

Storage instructionsStore at -20°C. Please refer to protocols.

Components	1000 tests
10X Solubilization Solution	1 x 22ml
20X Wash Buffer	1 x 50ml

Components	1000 tests
Doxorubicin	1 x 100µl
Fixation Solution II	1 x 55ml
SRB Solution	1 x 50ml

## Images



Dose-response curve of HepG2 (a), MCF7 (b) and HEK-293 (c) cells, after exposure to doxorubicin for 72 hours, as determined using the Sulforhodamine B Cell Cytotoxicity Assay Kit (Colorimetric). Assays were performed according to the kit protocol in triplicate.

**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