

Product datasheet

Cell Cycle Assay Kit (Fluorometric - Red) ab112117

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

Overview

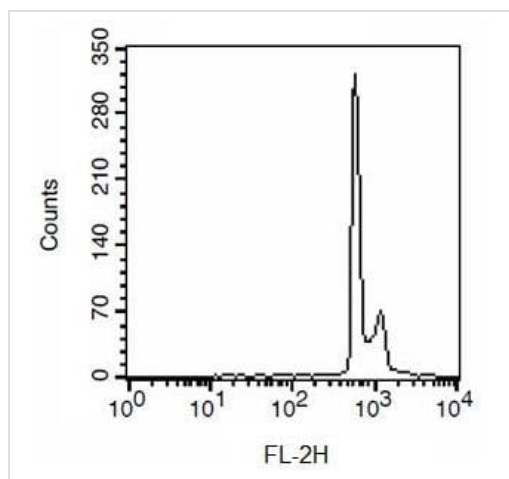
Product name	Cell Cycle Assay Kit (Fluorometric - Red)
Detection method	Fluorescent
Sample type	Adherent cells, Suspension cells
Assay type	Cell-based (quantitative)
Species reactivity	Reacts with: Mammals, Other species
Product overview	<p>ab112117 is designed to monitor cell cycle progression and proliferation by using our proprietary Nuclear Red CCS1 in fixed cells. The dye passes through a permeabilized membrane and intercalates into cellular DNA. The signal intensity of Nuclear Red CCS1 is directly proportional to DNA content after RNA is degraded by RNase provided in the kit. The percentage of cells in a given sample that are in G0/G1, S and G2/M phases, as well as the cells in the sub-G1 phase prior to apoptosis can be monitored with a flow cytometer at Ex/Em = 490/620 nm (FL2 channel).</p> <p>Visit our FAQs page for tips and troubleshooting.</p>
Notes	Store ab112117 desiccated.
Platform	Flow cytometer

Properties

Storage instructions Store at -20°C. Please refer to protocols.

Components	100 tests
100X RNase A	1 x 250µl
Assay Buffer	1 x 50ml
Nuclear Red CCS1	1 x 250µl

Images



Flow Cytometry - Cell Cycle Assay Kit (Fluorometric
- Red) (ab112117)

DNA profile in growing Jurkat cells.

Jurkat cells were dye-loaded with Red Fluorescence Probe and RNase A for 30 minutes. The fluorescence intensity of Red Fluorescence Probe was measured with a flow cytometer, using the FL2 channel.

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