

BrdU Cell Proliferation ELISA Kit (colorimetric) ab126556

★★★★★ [2 Abreviews](#) [160 References](#) [7 Images](#)

Overview

| | |
|------------------|--|
| Product name | BrdU Cell Proliferation ELISA Kit (colorimetric) |
| Detection method | Colorimetric |
| Sample type | Adherent cells, Suspension cells |
| Assay type | Sandwich (qualitative) |
| Sensitivity | < 40 cells/well |
| Assay time | 2h 30m |
| Assay duration | Multiple steps standard assay |
| Product overview | <p>BrdU Cell proliferation ELISA kit (ab126556) is an indirect ELISA kit for the detection of BrdU incorporation into newly synthesized DNA of actively proliferating cells. It involves incorporation of BrdU into cells cultured in microtiter plates using the cell layer as the solid phase. During the final 2 to 24 hours of culture BrdU is added to wells of the microtiter plate. BrdU will be incorporated into the DNA of dividing cells. To enable antibody binding to the incorporated BrdU cells must be fixed, permeabilized and the DNA denatured. This is all done in one step by treatment with Fixing Solution. Detector anti-BrdU monoclonal antibody is pipetted into the wells and allowed to incubate for one hour, during which time it binds to any incorporated BrdU. Unbound antibody is washed away and horseradish peroxidase-conjugated goat anti-mouse antibody is added, which binds to the Detector Antibody.</p> |

The horseradish peroxidase catalyzes the conversion of the chromogenic substrate tetra-methylbenzidine (TMB) from a colorless solution to a blue solution (or yellow after the addition of stopping reagent), the intensity of which is proportional to the amount of incorporated BrdU in the cells. The colored reaction product is quantified using a spectrophotometer.

The resultant assay is sensitive, rapid, easy to perform and applicable to high sample throughput. In addition to evaluation of cell proliferation, information such as cell number, morphology and analysis of cellular antigens can be obtained from a single culture.

Store kit at +4°C immediately upon receipt except for the Prediluted Anti-BrdU Detector Antibody and Peroxidase Goat anti-mouse IgG (2000X) which must be stored at -20°C.

Notes

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances. It is the responsibility of our customers to check the necessity of application of REACH

Authorisation, and any other relevant authorisations, for their intended uses.

Platform

Microplate

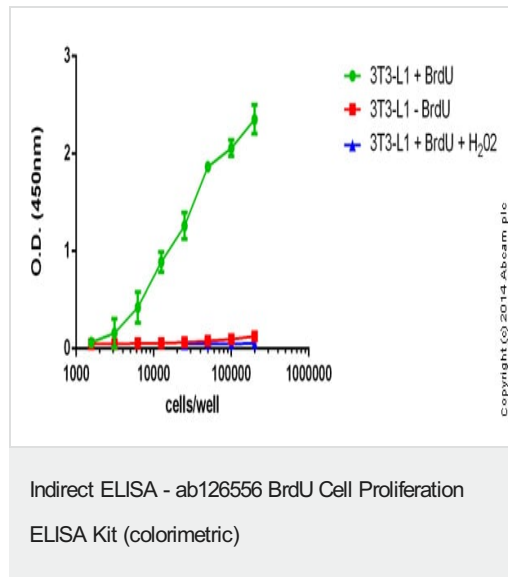
Properties

Storage instructions

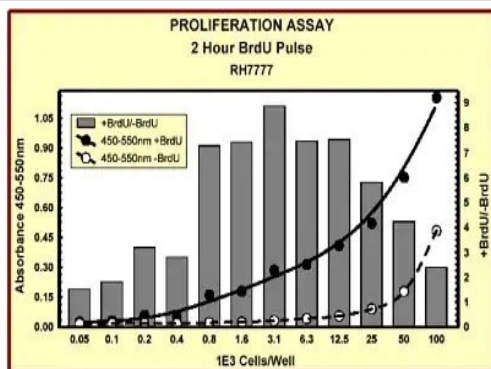
Store at +4°C. Please refer to protocols.

| Components | 200 tests |
|---|-----------|
| 500X BrdU Reagent | 1 x 15µl |
| Conjugate Diluent | 1 x 25ml |
| Fixing Solution | 2 x 20ml |
| Peroxidase Goat anti-mouse IgG (2000X) | 1 x 15µl |
| Plate wash concentrate (50X) | 1 x 90ml |
| Prediluted anti-BrdU detecting antibody | 1 x 20ml |
| Stop Solution | 1 x 25ml |
| TMB Substrate | 1 x 25ml |

Images

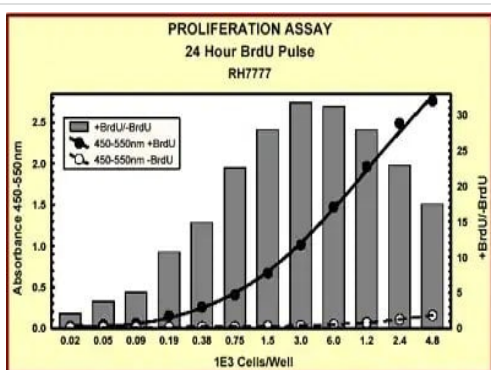


Cell proliferation measured in fibroblasts showing cell amounts vs. optical densities



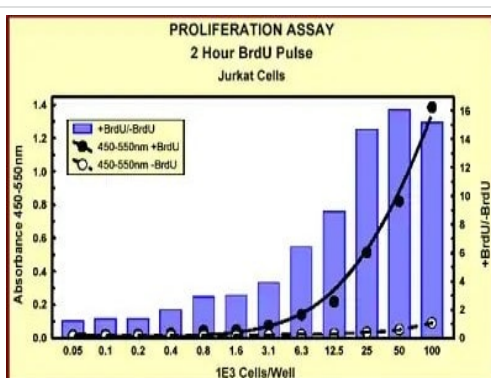
Indirect ELISA - BrdU Cell Proliferation ELISA Kit
(colorimetric) (ab126556)

Detection of RH7777 (adherent) cells per well with 2 hour pulse with BrdU. Y axis - left, OD 450-550 nm. Y axis-right, signal-to-noise ratio.



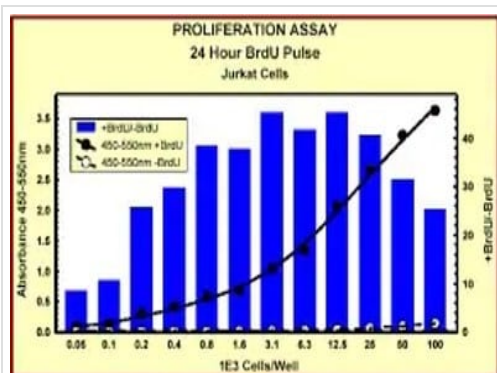
Indirect ELISA - BrdU Cell Proliferation ELISA Kit
(colorimetric) (ab126556)

Detection of RH7777 (adherent) cells per well with 24 hour pulse with BrdU. Y axis - left, OD 450-550 nm. Y axis-right, signal-to-noise ratio.



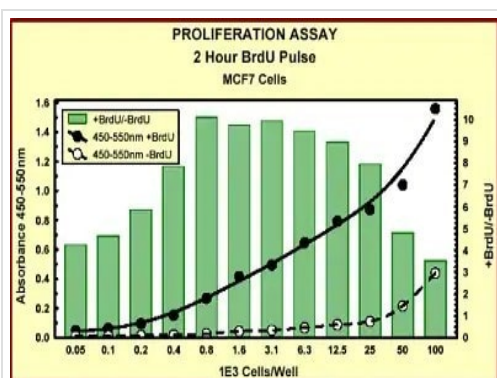
Indirect ELISA - BrdU Cell Proliferation ELISA Kit
(colorimetric) (ab126556)

Detection of Jurkat cells (non-adherent) per well with 2 hour pulse with BrdU. Y axis - left, OD 450-550 nm. Y axis-right, signal-to-noise ratio.



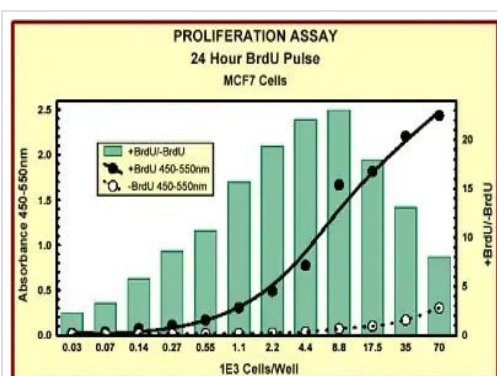
Indirect ELISA - BrdU Cell Proliferation ELISA Kit
(colorimetric) (ab126556)

Detection of Jurkat cells (non-adherent) per well with 24 hour pulse with BrdU. Y axis - left, OD 450-550 nm. Y axis-right, signal-to-noise ratio.



Indirect ELISA - BrdU Cell Proliferation ELISA Kit
(colorimetric) (ab126556)

Detection of MCF7 cells per well with 2 hour pulse with BrdU. Y axis - left, OD 450-550 nm. Y axis-right, signal-to-noise ratio.



Indirect ELISA - BrdU Cell Proliferation ELISA Kit
(colorimetric) (ab126556)

Detection of MCF7 cells per well with 24 hour pulse with BrdU. Y axis - left, OD 450-550 nm. Y axis-right, signal-to-noise ratio.

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