

BrdU Cell Proliferation ELISA Kit (chemiluminescent) ab126572

3 References

概述

产品名称	BrdU Cell Proliferation ELISA试剂盒(chemiluminescent)
检测方法	Luminescent
样品类型	Adherent cells, Suspension cells
检测类型	Sandwich (qualitative)
灵敏度	< 10 cells/well
检测时间	2h 00m
实验步骤	Multiple steps standard assay
产品概述	<p>ab126572 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> <p>The horseradish peroxidase catalyzes the oxidation of diacylhydrazides. The reaction product, in its excited state, decays yielding light in the process. This reaction has a large dynamic range spanning many logs of relative luminescence. The intensity of the luminescence is proportional to the amount of incorporated BrdU in the cells. The reaction is quantified using a luminometer and is plotted as relative light units per second (RLU/s).</p> <p>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.</p>
说明	<p>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.</p> <p>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.</p>
平台	Microplate

性能

存放说明

Please refer to protocols.

组件	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
Reaction Buffer	1 x 24ml
Substrate	1 x 800µl

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.cn/abpromise> or contact our technical team.

Terms and conditions

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