

Serine Protease Assay Kit (SR101-Leu-CMK) ab270776

医薬用外劇物

製品の概要

製品名	Serine Protease Assay Kit (SR101-Leu-CMK)
製品の概要	<p>These Serine Protease Assay Kits enable researchers to detect intracellular chymotrypsin-like serine protease activity <i>in vitro</i> without lysing the cell.</p> <p>FLISP® (Fluorescent Labeled Inhibitors of Serine Proteases) reagents are non-cytotoxic green or red fluorescent inhibitors that covalently bind with active serine protease enzymes. FLISP is a powerful method to assess the intracellular levels of chymotrypsin-like serine protease activity <i>in vitro</i>. Just add FLISP directly to the cell culture media, incubate, and wash. Because FLISP inhibitors are cell-permeant, they will efficiently diffuse in and out of all living cells. If there is an active chymotrypsin-like enzyme inside the cell, it will covalently bind with the FLISP inhibitor and retain the green or red fluorescent signal within the cell. Cells containing lower concentrations of chymotrypsin-like enzyme activity will retain a lower level of fluorescence compared to cells containing higher concentrations of this effector enzyme. There is no interference from pro-enzymes nor inactive forms of the enzymes. If the treatment is activating chymotrypsin-like serine proteases, positive cells will fluoresce brighter than the normal baseline negative cells, thus enabling researchers to clearly differentiate the populations.</p> <p>Produced using a non-baculovirus expression system.</p>
アプリケーション	適用あり: FM, Flow Cyt

製品の特性

保存方法 Please refer to protocols.

内容	25 tests	100 tests
10X Cellular Wash Buffer	1 x 15ml	1 x 60ml
Fixative	1 x 6ml	1 x 6ml
Hoechst 33342, 200 µg/mL	1 vial	1 vial
Red SR101-Leu-CMK-FLISP inhibitor reagent	1 vial	4 vials

アプリケーション

The Abpromise guarantee Abpromise保証は、次のテスト済みアプリケーションにおけるab270776の使用に適用されます
アプリケーションノートには、推奨の開始希釈率がありますが、適切な希釈率につきましてはご確認ください。

アプリケーション	Abreviews	特記事項
FM		Use at an assay dependent concentration.
Flow Cyt		Use at an assay dependent concentration.

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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