

Product datasheet

Serine Protease Assay Kit (FAM-Phe-DAP) ab270782

Overview

Product name	Serine Protease Assay Kit (FAM-Phe-DAP)
Product overview	<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>Fluorescent inhibitor reagents are non-cytotoxic green or red fluorescent inhibitors that covalently bind with active serine protease enzymes. This approach is a powerful method to assess the intracellular levels of chymotrypsin-like serine protease activity <i>in vitro</i>. Just add the fluorescent inhibitor reagent directly to the cell culture media, incubate, and wash. Because these 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 fluorescent inhibitor reagent 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>
Tested applications	Suitable for: Flow Cyt, FM

Properties

Storage instructions Please refer to protocols.

Components	25 tests	100 tests
10X Cellular Wash Buffer	1 x 15ml	1 x 60ml
Fixative	1 x 6ml	1 x 6ml
Green FAM-Phe-DAP fluorescent inhibitor reagent	1 vial	4 vials
Hoechst 33342, 200 µg/mL	1 vial	1 vial
Propidium Iodide, 250 µg/mL	1 x 1ml	1 x 1ml

Applications

The Abpromise guarantee Our [Abpromise guarantee](#) covers the use of ab270782 in the following tested applications. The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
Flow Cyt		Use at an assay dependent concentration.
FM		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.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