

## Datasheet: APO006

<b>Description:</b>	FLISP™ FAM-Phe-CMK SERINE PROTEASE ASSAY KIT
<b>Name:</b>	SERINE PROTEASE ASSAY KIT
<b>Format:</b>	FAM (Green Fluorescence)
<b>Product Type:</b>	Kits
<b>Quantity:</b>	25 TESTS

## Product Details

### Applications

This product has been reported to work in the following applications. This information is derived from testing within our laboratories, peer-reviewed publications or personal communications from the originators. Please refer to references indicated for further information. For general protocol recommendations, please visit [www.bio-rad-antibodies.com/protocols](http://www.bio-rad-antibodies.com/protocols).

	Yes	No	Not Determined	Suggested Dilution
Flow Cytometry	▪			
Immunofluorescence	▪			

Where this product has not been tested for use in a particular technique this does not necessarily exclude its use in such procedures. Suggested working dilutions are given as a guide only. It is recommended that the user titrates the product for use in their own system using appropriate negative/positive controls.

Max Ex/Em	Fluorophore	Excitation Max (nm)	Emission Max (nm)
	FAM	494	520

### Product Information

Fluorescent Labeled Inhibitors of Serine Protease (FLISP) kits allow the analysis of intracellular chymotrypsin-like serine protease activity in live cells. Serine protease activation plays major roles in several different cellular functions including apoptosis, tumor malignancy and as diagnostic and prognostic indicators of breast, head and neck carcinomas. Serine proteases are also thought to play a role in infections and transplant rejection.

### Test Principle

Active chymotrypsin-like enzymes within the cell bind covalently to the FLISP FAM-Phe-CMK inhibitor probe, retaining the green fluorescent signal within the cell. The higher the concentration of active chymotrypsin-like enzymes within the cell, the higher the level of fluorescence observed. FLISP kits are suitable for use with suspension or adherent cells and can be used with a flow cytometer, fluorescence microscope or a fluorescence plate reader.

FLISP inhibitors are cell-permeable and non-cytotoxic fluorochrome-labeled analogs of the first serine protease inhibitor, tosyl-phenylalanylchloromethyl ketone (TPCK), labeled with carboxyfluorescein (FAM) which fluoresces green. They are available with either a chloromethyl ketone (CMK) or diphenyl 1-(N-peptidylamino) alkanephosphonate (DAP) reactive group containing compound. Specificity for chymotrypsin-like enzymes is determined by the amino acid targeting peptide, the FAM-Phe-CMK probe inhibits chymotrypsin.

### Reagents In The Kit

FAM-Phe-CMK FLISP Reagent - lyophilized  
 10X Wash Buffer, 15 mL  
 Fixative, 6 mL

Propidium Iodide, 1 mL  
Hoechst Stain, 1 mL

---

<b>Instructions For Use</b>	Instructions for use can be found at <a href="http://www.bio-rad-antibodies.com/uploads/IFU/APO006.pdf">www.bio-rad-antibodies.com/uploads/IFU/APO006.pdf</a>
<b>Storage</b>	MULTIPLE STORAGE CONDITIONS APPLY ON ARRIVAL. Store the unopened kit (and each unopened component) according to the storage instructions on each component label. Store the FLISP reagent at -20°C. Once reconstituted in DMSO, the FLISP stock should be stored at -20°C for 6 months, protected from light. Avoid repeated freezing and thawing.
<b>Shelf Life</b>	Please see label for expiry date
<b>Acknowledgements</b>	FLISP is a trademark of Immunochemistry Technologies, LLC.
<b>Health And Safety Information</b>	Material Safety Datasheet documentation #20399 #10471 #10498 #10476 #10477 available at: FAM-Phe-CMK FLISP Reagent (20399): <a href="https://www.bio-rad-antibodies.com/uploads/MSDS/20399.pdf">https://www.bio-rad-antibodies.com/uploads/MSDS/20399.pdf</a> 10X Wash Buffer (10471): <a href="https://www.bio-rad-antibodies.com/uploads/MSDS/10471.pdf">https://www.bio-rad-antibodies.com/uploads/MSDS/10471.pdf</a> Fixative (10498): <a href="https://www.bio-rad-antibodies.com/uploads/MSDS/10498.pdf">https://www.bio-rad-antibodies.com/uploads/MSDS/10498.pdf</a> Hoechst Stain (10476): <a href="https://www.bio-rad-antibodies.com/uploads/MSDS/10476.pdf">https://www.bio-rad-antibodies.com/uploads/MSDS/10476.pdf</a> Propidium Iodide Reagent (10477): <a href="https://www.bio-rad-antibodies.com/uploads/MSDS/10477.pdf">https://www.bio-rad-antibodies.com/uploads/MSDS/10477.pdf</a>
<b>Regulatory</b>	For research purposes only

---

## Related Products

### Recommended Useful Reagents

[SR FLICA™ POLY CASPASE KIT \(ICT916\)](#)  
[SR FLICA™ POLY CASPASE KIT \(ICT917\)](#)  
[FAM FLICA™ POLY CASPASE KIT \(ICT091\)](#)  
[FAM FLICA™ POLY CASPASE KIT \(ICT092\)](#)  
[FLICA™ 660 POLY CASPASE KIT \(ICT9120\)](#)

<b>North &amp; South America</b>	Tel: +1 800 265 7376 Fax: +1 919 878 3751 Email: <a href="mailto:antibody_sales_us@bio-rad.com">antibody_sales_us@bio-rad.com</a>	<b>Worldwide</b>	Tel: +44 (0)1865 852 700 Fax: +44 (0)1865 852 739 Email: <a href="mailto:antibody_sales_uk@bio-rad.com">antibody_sales_uk@bio-rad.com</a>	<b>Europe</b>	Tel: +49 (0) 89 8090 95 21 Fax: +49 (0) 89 8090 95 50 Email: <a href="mailto:antibody_sales_de@bio-rad.com">antibody_sales_de@bio-rad.com</a>
----------------------------------	---	------------------	---	---------------	---

'M312675:180226'

Printed on 02 May 2018

---

© 2018 Bio-Rad Laboratories Inc | [Legal](#) | [Imprint](#)