

Datasheet: ICT096

Description:	FAM FLICA™ CASPASE-6 KIT
Name:	CASPASE-6
Format:	FAM (Green Fluorescence)
Product Type:	Kits
Quantity:	100 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.

	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.

Product Information

The FAM FLICA™ Caspase 6 kit uses a quick and easy method to analyze active caspases in apoptotic cells.

Test Principle

Caspase FLICA™ kits measure apoptosis by detecting active caspases in whole, living cells. These kits do not work by using antibodies or as an ELISA. Instead, their methodology is based on a unique cell-permeable and non-cytotoxic reagent called the Fluorochrome Inhibitor of Caspases (FLICA). The FLICA™ reagent contains a caspase inhibitor sequence linked to a green (Carboxyfluorescein, FAM) fluorescent probe.

The Caspase FLICA™ Kits are suitable for cells in suspension, adherent cells, thin tissue sections (but not fixed or paraffin-embedded cells) from many species including mammalian, insect and yeast. Different cell types, e.g. HeLa, primary neurons, macrophages and lymphocytes have also been successfully studied with these kits.

This kit can be used with a flow cytometer, fluorescence microscope or a fluorescence plate reader using black microtitre plates.

Reagents In The Kit

4 vials of FAM-VEID-FMK FLICA™ Reagent - lyophilized
10X Wash Buffer, 60 mL
Fixative, 6 mL
Propidium Iodide, 1 mL
Hoechst Stain, 1 mL

Instructions For Use

Instructions for use can be found at www.bio-rad-antibodies.com/uploads/IFU/ICT096.pdf

Storage	Store the unopened kit (and each unopened component) at +4°C until the expiration date. Protect the FLICA™ reagent from light at all times. Once reconstituted, the 150X FLICA™ stock should be stored at -20°C protected from light.
Shelf Life	Please see label for expiry date.
Acknowledgements	FLICA™ is a trademark of Immunochemistry Technologies, LLC.
Health And Safety Information	Material Safety Datasheet documentation #20283 #10471 #10498 #10476 #10477 available at: FAM-VEID-FMK FLICA²/sup> Reagent (20283): https://www.bio-rad-antibodies.com/uploads/MSDS/20283.pdf 10X Wash Buffer (10471): https://www.bio-rad-antibodies.com/uploads/MSDS/10471.pdf Fixative (10498): https://www.bio-rad-antibodies.com/uploads/MSDS/10498.pdf Hoechst Stain (10476): https://www.bio-rad-antibodies.com/uploads/MSDS/10476.pdf Propidium Iodide Reagent (10477): https://www.bio-rad-antibodies.com/uploads/MSDS/10477.pdf
Regulatory	For research purposes only

North & South America	Tel: +1 800 265 7376 Fax: +1 919 878 3751 Email: antibody_sales_us@bio-rad.com	Worldwide	Tel: +44 (0)1865 852 700 Fax: +44 (0)1865 852 739 Email: antibody_sales_uk@bio-rad.com	Europe	Tel: +49 (0) 89 8090 95 21 Fax: +49 (0) 89 8090 95 50 Email: antibody_sales_de@bio-rad.com
----------------------------------	---	------------------	---	---------------	---

'M310050:170914'

Printed on 01 May 2018