

## Datasheet: ICT095

|                      |                          |
|----------------------|--------------------------|
| <b>Description:</b>  | FAM FLICA™ CASPASE-6 KIT |
| <b>Name:</b>         | CASPASE-6                |
| <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.

### 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/ICT095.pdf](http://www.bio-rad-antibodies.com/uploads/IFU/ICT095.pdf)

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

## Related Products

### Recommended Useful Reagents

[FAM FLICA™ CASPASE-6 KIT \(ICT096\)](#)

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

'M310049:170914'

Printed on 01 May 2018