

Datasheet: ICT935

Description:	MAGIC RED™ CASPASE-3/7 KIT
Name:	CASPASE-3/7
Format:	MAGIC RED™ (Red Fluorescence)
Product Type:	Kits
Quantity:	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.

	Yes	No	Not Determined	Suggested Dilution
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 Magic Red™ Caspase 3 & 7 kit uses a quick and easy method to analyze intracellular active caspases in apoptotic cells.

Test Principle

Caspase Magic Red™ 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 cell-permeable and non-cytotoxic reagent which is cleaved in the presence of caspases to produce a fluorescent product. The Magic Red™ reagent contains a caspase inhibitor sequence (DEVD)₂ linked to a red (Cresyl Violet) fluorescent probe.

The Caspase Magic Red™ kits are suitable for cells in suspension, adherent cells, thin frozen tissue sections (but not fixed or paraffin-embedded cells) from many species including mammalian, insect and yeast. Different cell types, e.g. Jurkat, HL-60, THP-1, fibroblasts, UMUC-3, MCF-7 and U937 cells have also been successfully studied with these kits.

This kit can be used in fluorescence microscopy or with fluorescence plate readers using black microtitre plates.

Reagents In The Kit

Magic Red™ MR-(DEVD)₂ Reagent - lyophilized
Hoechst 33342 Stain, 1 mL
Acridine orange (AO), 0.5 mL

Instructions For Use

Instructions for use can be found at www.bio-rad-antibodies.com/uploads/IFU/ICT935.pdf

References

1. Wei, Z. *et al.* (2011) Notch activation enhances the microglia-mediated inflammatory response associated with focal cerebral ischemia. [Stroke](#). 42 (9): 2589-94.
2. Shibata, Y. *et al.* (2016) Controlled in-cell release of caspase from photodegradable

nanoparticles using the PARCEL method [Applied Spectroscopy Reviews](#). Mar 13 [Epub ahead of print]

Storage	Store the unopened kit (and each unopened component) at +4°C until the expiration date. Protect the MR-(DEVD) ₂ reagent from light at all times. Once reconstituted, the 155X MR-(DEVD) ₂ stock should be stored at -20°C protected from light.
Shelf Life	Please see label for expiry date.
Acknowledgements	Magic Red™ is a trademark of Immunochemistry Technologies, LLC.
Health And Safety Information	Material Safety Datasheet documentation #20262 #10476 #10478 available at: Magic Red [®] /sup> MR-(DEVD) ₂ Reagent (20262): https://www.bio-rad-antibodies.com/uploads/MSDS/20262.pdf Hoechst Stain (10476): https://www.bio-rad-antibodies.com/uploads/MSDS/10476.pdf Acridine Orange Stain (10478): https://www.bio-rad-antibodies.com/uploads/MSDS/10478.pdf
Regulatory	For research purposes only

Related Products

Recommended Useful Reagents

[MAGIC RED™ CASPASE-3/7 KIT \(ICT936\)](#)

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
----------------------------------	---	------------------	---	---------------	---

'M289540:160616'

Printed on 30 Apr 2018