

## Datasheet: ICT936

<b>Description:</b>	MAGIC RED™ CASPASE-3/7 KIT
<b>Name:</b>	CASPASE-3/7
<b>Format:</b>	MAGIC RED™ (Red Fluorescence)
<b>Product Type:</b>	Kits
<b>Quantity:</b>	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](http://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)<sub>2</sub> 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)<sub>2</sub> 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/ICT936.pdf](http://www.bio-rad-antibodies.com/uploads/IFU/ICT936.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 \[Epub ahead of print\]](#)

---

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

---

**North & South** Tel: +1 800 265 7376

**America** Fax: +1 919 878 3751

Email: [antibody\\_sales\\_us@bio-rad.com](mailto: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](mailto: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](mailto:antibody_sales_de@bio-rad.com)

'M289541:160616'

**Printed on 30 Apr 2018**

---

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