

## Product datasheet

# Autophagy Assay Kit (Red) ab270790

### Overview

**Product name** Autophagy Assay Kit (Red)

**Product overview** Autophagy Assay kit enables researchers to detect and monitor the in vitro development of autophagy in living cells. Autophagy Probe, Red is a cell-permeant aliphatic molecule that fluoresces brightly when inserted in the lipid membranes of autophagosomes and autolysosomes. Autophagy Probe, Red can be readily detected by flow cytometry with optimal excitation at 590 nm and peak emission at 620 nm. Autophagy Probe, Red is for research use only. Not for use in diagnostic procedures.

Produced using a non-baculovirus expression system.

### Properties

**Storage instructions** Please refer to protocols.

Components	50 tests	200 tests
10X Cellular Assay Buffer	1 x 15ml	1 x 60ml
Autophagy Probe, Red, 50 tests per vial	1 vial	4 vials
Fixative	1 x 6ml	1 x 6ml

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

### Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

## **Terms and conditions**

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors