

Version 1 Last updated 30 January 2020

ab270786

Caspase 8 Assay Kit (Red)

View ab270786

Caspase 8 Assay Kit (Red) datasheet:

www.abcam.com/ab270786

(use www.abcam.cn/ab270786 for China, or www.abcam.co.jp/ab270786 for Japan)

For the detection of caspase activity in cultured cells and tissues.

This product is for research use only and is not intended for diagnostic use.

Table of Contents

1. Overview	1
2. Materials Supplied and Storage	2
3. Materials Required, Not Supplied	3
4. General guidelines, precautions, and troubleshooting	4
5. Experimental Preparation	5
6. Controls	6
7. Apoptosis Induction	7
8. Preparation of SR-LETD-FMK	7
9. Preparation of 1X Apoptosis Wash Buffer	8
10. Hoechst 33342	8
11. Fixative	9
12. Staining Protocol for Suspension Cells	9
13. Staining Protocol for Adherent Cells	11
14. Microscopy analysis	13
15. Fluorescent Plate Reader Analysis	14
16. Flow Cytometry Analysis	15
17. Notes	17

1. Overview

Caspases play important roles in apoptosis and inflammation. Caspase 8 Assay Kit (Red) (ab270786) is used by researchers seeking to quantitate apoptosis via caspase activity in cultured cells and tissues.

The reagent SR-LETD-FMK enters each cell and irreversibly binds to activated caspase-8. Because the SR-LETD-FMK reagent becomes covalently coupled to the active enzymes, it is retained within the cell, while any unbound SR-LETD-FMK reagent diffuses out of the cell and is washed away. The remaining red fluorescent signal is a direct measure of the active caspase-8 enzyme activity present in the cell at the time the reagent was added. Cells that contain the bound SR-LETD-FMK can be analyzed by a fluorescence plate reader, fluorescence microscopy, or flow cytometry. Cells labeled with the SR-LETD-FMK reagent may be read immediately or preserved for 16 hours using the fixative included in the kit. Unfixed samples may also be analyzed with Hoechst 33342 to detect changes in nuclear morphology.

There is no interference from pro-caspases or inactive forms of the enzymes. If the treatment is causing cell death via apoptosis, apoptotic cells will have an elevated level of caspase activity relative to nonapoptotic or negative control cells and fluoresce with SR-LETD-FMK.

2. Materials Supplied and Storage

Store the kit at +4°C upon receipt.

Δ Note: Once reconstituted with DMSO, use SR-LETD-FMK immediately, or store at $\leq -20^{\circ}\text{C}$ for up to 6 months, protected from light. Avoid repeated freeze/thaw cycles.

Item	Quantity	Storage Temperature
10X Apoptosis Wash Buffer (25 tests)	1 x 15 mL	+4°C
10X Apoptosis Wash Buffer (100 tests)	1 x 60 mL	
Fixative	1 x 6 mL	+4°C
Caspase inhibitor SR-LETD-FMK (25 tests)	1 vial	+4°C
Caspase inhibitor SR-LETD-FMK (100 tests)	4 vials	+4°C
Hoechst 33342	1 vial	+4°C

3. Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully perform this assay:

- DMSO, 50 μ L per vial to reconstitute SR-LETD-FMK
- DiH_2O , 135-540 mL to dilute 10X Apoptosis Wash Buffer
- Phosphate buffered saline (PBS) pH 7.4, up to 100 mL, to dilute SR-LETD-FMK and handle cells
- FBS and/or BSA to add to the buffer when handling cells
- Cultured cells treated with the experimental conditions ready to be labeled
- Reagents to induce apoptosis or trigger caspase activity, such as staurosporine or camptothecin
- Hemocytometer
- Centrifuge at 200 $\times g$
- 15 mL polystyrene centrifuge tubes (1 per sample)
- Fluorescence microscope
- Flow Cytometer
- Fluorescent plate reader

4. General guidelines, precautions, and troubleshooting

Please observe safe laboratory practice and consult the safety datasheet.

For general guidelines, precautions, limitations on the use of our assay kits and general assay troubleshooting tips, particularly for first time users, please consult our guide:

www.abcam.com/assaykitguidelines

For typical data produced using the assay, please see the assay kit datasheet on our website.

5. Experimental Preparation

- 5.1 Staining apoptotic cells with SR-LETD-FMK can be completed within a few hours. However, SR-LETD-FMK is used with living cells, which require periodic maintenance and cultivation several days in advance.
- 5.2 In addition, once the proper number of cells has been cultivated, time must be allotted for the experimental treatment or apoptosis induction process which typically requires a 2-6 hour incubation at 37°C based on the cell line and concentration. Create cell populations, such as:
 - a. Cells that were exposed to the experimental condition or treatment
 - b. A placebo population of cells that received a blank treatment instead of the experimental treatment.
- 5.3 As SR-LETD-FMK detects the presence of catalytically active forms of caspase enzymes, plan the experiment so that SR-LETD-FMK will be diluted and administered at the time when caspases are expected to be activated in the cells.
- 5.4 The recommended volume of 30-60X SR-LETD-FMK is 5-10 μL per 300 μL of cells at $3-5 \times 10^5$ cells/mL, but the amount may vary based on the experimental conditions and the instrument used for analysis. Each investigator should adjust the amount of SR-LETD-FMK to accommodate the particular cell line and research conditions.
- 5.5 Culture cells to a density optimal for the specific experiment or apoptosis induction protocol. Cell density should not exceed 10^6 cells/mL. Cells cultivated in excess of this concentration may begin to naturally enter apoptosis.
- 5.6 An initial experiment may be necessary to determine when and how much SR-LETD-FMK to use as the resulting positive signal is a direct measurement of caspase activity occurring during the incubation period.

6. Controls

Create experimental samples and control cell populations:

- Treated experimental population(s): cells exposed to the experimental condition(s).
- Negative control: non-treated cells grown in a normal cell culture environment.
- Positive control: cells induced to undergo apoptosis using a known apoptosis induction protocol.

The induced positive cell population and negative control cell population tubes should come from a common pool of cells and contain similar quantities of cells. Create negative controls by culturing an equal volume of noninduced cells for every labeling condition. For example, if labeling with SR-LETD-FMK and Hoechst 33342 (which is optional), make eight populations:

Control #	
1 & 2	Unlabeled: induced and non-induced
3 & 4	SR-LETD-FMK -labeled: induced and non-induced
5 & 6	SR-LETD-FMK -labeled and Hoechst-labeled: induced and non-induced
7 & 8	Hoechst-labeled: induced and non-induced

7. Apoptosis Induction

Prior to commencing the experiment, determine a reproducible method for obtaining a positive control by triggering caspase activity. This process varies significantly with each cell line. For example, apoptosis may be induced with 2-4 $\mu\text{g}/\text{mL}$ camptothecin or 1-2 μM staurosporine for >4 hours.

8. Preparation of SR-LETD-FMK

SR-LETD-FMK is supplied as a lyophilized powder that may be slightly visible as an iridescent sheen inside the vial. Protect from light and use gloves when handling. **Because the 30-60X SR-LETD-FMK solution must be used immediately, prepare it just before staining.**

- 8.1 Reconstitute each vial of SR-LETD-FMK with 50 μL DMSO to form the 150-300X stock concentrate. The stock solution should be pink to red in color. Once reconstituted, it may be stored at $\leq -20^{\circ}\text{C}$ for 6 months protected from light and thawed no more than twice during that time.
- 8.2 Immediately prior to addition to the samples and controls, dilute SR-LETD-FMK 1/5 by adding 200 μL PBS to each vial to form the 30-60X SR-LETD-FMK solution. **Use 30-60X SR-LETD-FMK within 30 minutes of dilution into aqueous buffers.**

9. Preparation of 1X Apoptosis Wash Buffer

10X Apoptosis Wash Buffer is an isotonic solution used to wash cells following exposure to SR-LETD-FMK. It contains mammalian proteins to stabilize cells stained with SR-LETD-FMK and sodium azide to retard contamination (1X Apoptosis Wash Buffer contains 0.01% w/v sodium azide). Cell media containing FBS and other additives may be used instead of 1X Apoptosis Wash Buffer.

- 10X Apoptosis Wash Buffer may form precipitates during cold storage. If this happens, gently warm it until all crystals have dissolved. Do not boil.
- Dilute 10X Apoptosis Wash Buffer 1/10 in diH₂O. For example, add 15 mL 10X Apoptosis Wash Buffer to 135 mL diH₂O for a total of 150 mL.

Δ Note: 1X Apoptosis Wash Buffer may be stored at 4°C and used within 1 week or frozen and used within 6 months.

10. Hoechst 33342

Hoechst 33342 is a cell-permeant nuclear stain that emits blue fluorescence when bound to double stranded DNA. It is used to stain the nuclei of living or fixed cells, to distinguish condensed pyknotic nuclei in apoptotic cells, and for cell cycle studies.

Hoechst 33342 is provided ready-to-use at 200 µg/mL. Hoechst 33342 can be used with SR-LETD-FMK to label the nuclei of live, dying, and apoptotic cells.

When bound to nucleic acids, it has a maximum absorbance at 350 nm and a maximum emission at 480 nm. It is revealed under a microscope using a UV-filter with excitation at 365 nm and emission at 480 nm.

Δ Note: Hoechst 33342 contains a low concentration of Bis benzimide H 33342 trihydrochloride (CAS 23491-52-3) which is below the threshold for reporting on the safety data sheet (SDS). It is a suspected mutagen at high concentrations. Prolonged skin contact may cause redness and irritation. Because of the small quantity of product, the health hazard is small. See SDS for further information.

11. Fixative

The Fixative is a formaldehyde solution designed to cross-link and aggregate intracellular components. If the stained cell populations cannot be evaluated immediately after labeling with SR-LETD-FMK, add Fixative at a ratio of 1:5-1:10. For example, to use Fixative at 1:10, add 100 μ L Fixative to 900 μ L cells. Never add Fixative until all the staining and final wash steps have been completed. Fixed cells may be stored protected from light on ice or at 2-8°C for up to 16 hours.

The fixative will not interfere with the sulforhodamine (SR) label. Do not use absolute ethanol- or methanol-based fixatives, as they may inactivate the SR-LETD-FMK label.

Δ Note: Fixative contains formaldehyde <10% and methanol <5% and is harmful. Avoid contact with skin, eyes and clothing by wearing lab coat, gloves and safety glasses. In case of exposure, immediately flush eyes or skin with water. See SDS for further information.

12. Staining Protocol for Suspension Cells

Prepare experimental and control cell populations. Ideally, cell concentration should be 3-5 x 10⁵ cells/mL. The concentration should not exceed 10⁶ cells/mL, as cells cultivated in excess of this concentration may begin to naturally enter apoptosis. Just prior to staining with SR-LETD-FMK, cells may need to be concentrated to 2-5 x 10⁶ cells/mL as both microscopy and plate reader analysis methods (Sections 14 and 15) require high cell concentrations. Start with a larger volume of cells at 3-5 x 10⁵ cells/mL (which is a typical density for cell culture) and then concentrate cells and resuspend to 300 μ L per sample when ready for SR-LETD-FMK staining.

- 12.1 Expose cells to the experimental and control conditions. Include the following controls:
 - a. Positive control cells induced to undergo apoptosis.
 - b. Negative control cells not induced to undergo apoptosis.
- 12.2 If analyzing with a fluorescence microscope or plate reader, concentrate cells to 2-5 x 10⁶ cells/mL just prior to SR-LETD-FMK staining. Fluorescence microscopy requires an excess of 2 x 10⁶

cells/mL to obtain 5-20 cells per image field. Flow cytometry can efficiently analyze samples at $3\text{-}5 \times 10^5$ cells/mL.

- 12.3** Transfer 290-295 μL cells into fresh tubes.
- 12.4** Add 5-10 μL 30-60X SR-LETD-FMK solution, forming a final volume of 300 μL . If different cell volumes were used, add SR-LETD-FMK at a ratio of 1:30-1:60. Mix by gently flicking the tubes. The amount of SR-LETD-FMK should be optimized for each cell line and experimental condition. However, 1:60 is generally sufficient for flow cytometry applications, and 1:30 is recommended for fluorescence microscopy or plate reader applications.
- 12.5** Incubate cells at 37°C protected from light. The incubation period may range from 30 minutes to several hours and should be optimized for each cell line and experimental condition. As cells may settle on the bottom of the tubes, gently resuspend by swirling cells every 10-20 minutes to ensure an even distribution of SR-LETD-FMK.
- 12.6** If cells are to be analyzed with a microscope, cells may be dually stained with Hoechst. Do not stain with Hoechst if using a plate reader (Section 15) or flow cytometer (Section 16). Add Hoechst at 0.5% v/v and incubate 5 minutes at 37°C . For example, if the cell suspension is at 300 μL , add 1.5 μL Hoechst.
- 12.7** Add 2 mL 1X Apoptosis Wash Buffer and gently mix.
- 12.8** Centrifuge at $200 \times g$ for 5-10 minutes at room temperature (RT).
- 12.9** Carefully remove and discard supernatants. Gently vortex pellets to disrupt clumping. Resuspend in 1 mL 1X Apoptosis Wash Buffer and gently mix.
- 12.10** Centrifuge cells at $200 \times g$ for 5-10 minutes at RT.
- 12.11** Carefully remove and discard supernatants. If analyzing by fluorescence microscopy or fluorescence plate reader, repeat wash process a third time. If using a flow cytometer, two wash steps are generally sufficient.
- 12.12** Gently vortex pellets to disrupt clumping.
 - If analyzing with a fluorescence microscope, go to Section 14.
 - If using a fluorescence plate reader, go to Section 15.
 - If using a flow cytometer, go to Section 19.

13. Staining Protocol for Adherent Cells

Adherent cells need to be handled carefully to avoid the loss of any cells that round up and come off the culture surface. In microscopy or plate reader applications where trypsinization is not required, adherent cells can be stained and washed directly on the chamber slide, well, or culture surface. To avoid losing cells that are no longer adherent during washing, spin down all overlay media and wash buffer and recombine washed cell pellets with the adherent samples prior to analysis.

Cells may be trypsinized to create suspensions, which may be labeled with SR-LETD-FMK before or after trypsinization. Avoid trypsinizing cells prior to labeling with a live/dead DNA dye, like 7-AAD. Cell membranes exposed to trypsin could be transiently permeable to live/dead dyes for a variable time depending upon the cell line.

- 13.1** Expose cells to the experimental and control conditions. Include the following controls:
- 13.2** If staining cells while adherent, go to Step 13.4. If suspension cells are required for the final analysis:
 - Remove overlay media. Spin to pellet any loose cells.
 - Trypsinize adherent cells. Alternatively, SR-LETD-FMK labeling can be performed first, followed by washing and trypsinization steps.
 - Neutralize with trypsin inhibitor, as found in cell culture media with 20% FBS.
 - Add 2-5 mL media
- 13.3** Prepare trypsinized cells for staining:
 - Centrifuge at 200 x g for 5-10 minutes at RT.
 - Remove all but ~100 μ L supernatant.
 - Resuspend cells in 300-500 μ L cell culture media containing 10-20% FBS.
 - If necessary, count cells and adjust the concentration and volume of cell suspension to fit the experiment
- 13.4** Add the SR-LETD-FMK working solution to samples at a v/v ratio of 1:30-1:60 and mix the cell suspension to disperse the SR-LETD-FMK reagent. If staining adherent cells, add SR-LETD-FMK directly to the overlay media. The concentration of SR-LETD-

FMK should be optimized for each cell line, experimental condition, and method of analysis:

- Flow cytometry analysis may provide the sensitivity to detect SR-LETD-FMK when used at 1:60.
 - For analysis by fluorescence microscopy, fluorescence plate reader, or applications where a higher staining concentration is needed, it is recommended that the SR-LETD-FMK working solution be used at 1:30. For example, if staining cells at 1:30, add 10 μ L SR-SR-LETD-FMK working solution to 290 μ L cells, forming a final volume of 300 μ L.
- 13.5** Incubate 30-60 minutes at 37°C, mixing gently every 10-20 minutes to disperse the reagent.
- 13.6** Wash cells. For trypsinized adherent (suspension) cells, go to Step 13.7. For adherent cells, go to Step 13.8.
- 13.7** Wash trypsinized adherent (suspension) cells:
- Add 2 mL 1X Apoptosis Wash Buffer and gently mix.
 - Centrifuge at 200 x *g* for 5-10 minutes at RT.
 - Carefully aspirate supernatant.
 - Resuspend samples in 1-2 mL wash buffer and gently mix.
 - Centrifuge a second time at 200 x *g* for 5-10 minutes at RT.
 - Carefully aspirate supernatant.
 - For flow cytometry analysis, two wash steps are generally sufficient. For microscopy or fluorescence plate reader analysis, repeat wash procedure a third time (resuspend samples, gently pellet by centrifugation, and carefully remove supernatants). Go to Step 13.9.
- 13.8** Wash adherent cells. To avoid losing cells that are no longer adherent during washing, spin down all overlay media and wash buffer and recombine washed cell pellets with the adherent samples prior to analysis.
- Carefully remove overlay media containing SR-LETD-FMK and replace with 1X Apoptosis Wash Buffer.
 - Incubate 10 minutes at 37°C to allow any unbound SR-LETD-FMK to diffuse out of cells.
 - Carefully remove and replace wash buffer with fresh wash buffer and incubate another 10 minutes at 37°C.
 - Gently remove overlay buffer and replace for a third wash step. Incubate 10 minutes at 37°C.

- Gently remove overlay buffer. Go to Step 13.9
- 13.9** If desired and using a microscope, cells may be counter-stained with ancillary dyes like the nuclear stain Hoechst 33342 or other compatible fluorescent dye.
- Resuspend cells or replace overlay buffer, add Hoechst 33342 at 0.5% v/v, and incubate 5 minutes at 37°C. For example, if the cell suspension or overlay volume is at 1 mL, add 5 µL Hoechst 33342.
 - Live/dead cell stains should not be used after trypsinization. Cell membranes exposed to trypsin could be transiently permeable to live/dead dyes for a variable time depending upon the cell line. To identify dead cells with a live/dead stain, perform the SR-LETD-FMK and ancillary dye staining and wash steps prior to using trypsin. To avoid false positives, include another wash step to remove excess live/dead dye prior to trypsinization.
 - If using any compatible ancillary dyes, follow the manufacturer's specific instructions for staining samples.
- 13.10** Read cells within 4 hours or fix.
- If analyzing with a fluorescence microscope, go to Section 14.
 - If using a fluorescence plate reader, go to Section 15.
 - If using a flow cytometer, go to Section 16.

14. Microscopy analysis

Follow Section 12 or Section 13.

- 14.1** Resuspend cells or replace overlay media in 300-500 µL 1X Apoptosis Wash Buffer and place on ice. At this point, the cells may be stained with other dyes, fixed for future viewing (Step 14.2), or observed immediately (Step 14.3).
- 14.2** If not viewing immediately, cells may be fixed for viewing up to 16 hours later.
- Add Fixative at a v/v ratio of 1:5-1:10.
 - Incubate 15 minutes at RT in the dark.
 - Place cells on a microscope slide and allow to dry.

- Briefly rinse cells with PBS.
 - Cover with mounting media and coverslip.
 - Store slides at 4°C for up to 16 hours
- 14.3** To view cells immediately, place 1 drop of cell suspension onto a microscope slide and cover with a coverslip.
- 14.4** Observe cells with a fluorescence microscope using a bandpass filter (excitation 550 nm, emission >580 nm) to view red fluorescence. Cells bearing active caspase enzymes covalently coupled to SR-LETD-FMK appear red. Hoechst 33342 can be seen using a UV-filter with excitation at 365 nm and emission at 480 nm.

15. Fluorescent Plate Reader Analysis

Follow Section 12 or Section 13, but omit optional nuclear staining with Hoechst 33342.

- 15.1** Determine the concentration and compare the cell density of each sample. The non-induced population may have more cells than the induced population, as some apoptotic cells in the induced samples may be lost during the wash steps. Adjust the volume of the cell suspensions to equalize the cell density. When ready to read, cells should be $>3 \times 10^6$ cells/mL. Adherent cells should be cultured to 80-90% confluency. Please note that some cell lines will not tolerate confluency levels $>60\%$; adjust as necessary for the particular cells being used.
- 15.2** If using suspension cells, pipette 100 μ L stained and washed cells per well into a black microtiter plate. Do not use clear plates. If using a bottom-reading instrument, use a plate with black walls and a clear bottom. Analyze at least 2 aliquots per sample. Avoid bubbles.
- 15.3** Perform an endpoint read. Set the excitation wavelength at 550 nm and the emission wavelength at 595 nm; if possible, use a 570 nm cut-off filter. SR-LETD-FMK is excited at 550-580 nm (excitation peak is 570 nm) and the emission optima is at 590-600 nm (emission peak is 590 nm).

16. Flow Cytometry Analysis

Flow cytometry with SR-LETD-FMK is best performed using a laser that can efficiently excite the fluorophore, such as a yellow-green 561 nm laser. A green 532 nm laser with a 610/20 filter pairing has also been shown to be effective. If yellow-green or green laser options are unavailable, a standard 488 nm blue excitation laser paired with an appropriate emission filter set, such as 585/40 (often FL-2/PE channel), can be used.

When using SR-LETD-FMK in combination with other fluorochromes for bi-color analyses, compensation may be necessary. Correcting for fluorescence spill over is especially important when using a sub-optimal excitation laser to excite SR-LETD-FMK in combination with a fluorochrome with significant spectral overlap with the SR-LETD-FMK detector. For instance, some green dyes, like FITC or Green Live/Dead Stain (catalog #6342), can spill over from FL-1 into FL-2. If using such a dye to co-label SR-LETD-FMK stained cells, consider reading the SR-LETD-FMK reagent in an alternate channel with reduced spectral overlap, such as 670LP (often FL-3/PerCP), or similar.

Because of its orangy-red profile, SR-LETD-FMK can be measured in flow cytometry with the red live/dead stain, 7-AAD (catalog #6163), to detect necrosis concurrently (Figure 7). Careful compensation and gating of SR-LETD-FMK and 7-AAD fluorophores (using the FL-2 and FL-3 channels, respectively) will distinguish the red 7-AAD live/dead fluorescence signal from the orangy-red SR-LETD-FMK caspase positive signal within each sample. This will reveal four populations of cells:

- Live unstained cells display only normal background or autofluorescence (revealed in lower left quadrant).
- Early apoptotic cells fluoresce orangy-red with SR-LETD-FMK (revealed in lower right quadrant).
- Late apoptotic cells are dually stained with SR-LETD-FMK and 7-AAD: they fluoresce orangy-red (they have active caspases) and red (the cell membrane has permeabilized). This population will be revealed in the upper right quadrant.
- Necrotic cells fluoresce red with 7-AAD (revealed in the upper left quadrant).

Follow Section 12 or Section 13, but omit optional nuclear staining with Hoechst 33342

- 16.1** Resuspend cells in 300 μ L 1X Apoptosis Wash Buffer and place on ice.
- 16.2** Cells may be fixed for analysis up to 16 hours later. Add Fixative at a v/v ratio of 1:5-1:10. Store samples at 2-8°C and protected from light.
- 16.3** Run the unstained control. If possible, adjust voltages to place the unstained sample in the first decade of the FL dot plots.
- 16.4** For single-color analysis, a 488 nm blue argon laser or comparable can be used with the emission filter pairing that best approximates 585/40 (often FL-2/PE channel). If other options are available, use the laser/ filter pairing that most closely resembles the excitation and emission optima of sulforhodamine B, i.e. 570 nm and 590-600 nm, respectively.
- 16.5** Generate a histogram with the log FL-2 on the X-axis versus the number of cells on the Y-axis. Caspase negative (SR-LETD-FMK -) cells will fall within the lower log fluorescence output decades of the FL-2 X-axis, whereas caspase-positive (SR-LETD-FMK +) cells will appear as a shoulder or as a separate peak on the right side of the negative peak histogram.
- 16.6** For bi-color analyses, run each single color control. Adjust compensation to remove spectral overlap from interfering FL channels. When the data have been correctly compensated, the median fluorescence intensity (MFI) values in non-primary detectors of any given single-stained control sample should be the same as an unstained control sample (e.g. an SR-LETD-FMK stained sample being read in FL-2 should have the same MFI in FL-3 as an unstained sample). For example, if reading SR-LETD-FMK in FL-2 and 7-AAD in FL-3:
 - Subtract a percentage of the fluorescence in the SR-LETD-FMK channel from the fluorescent channel used for 7-AAD (e.g. FL-3 - %FL-2).
 - Subtract a percentage of the fluorescence in the channel used for 7-AAD from the fluorescence in the SR-LETD-FMK channel (e.g. FL-2 - %FL-3).
- 16.7** Run the bi-color experimental samples and analyze.

17. Notes

Technical Support

Copyright © 2020 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

Austria

wissenschaftlicherdienst@abcam.com | 019-288-259

France

supportscientifique@abcam.com | 01.46.94.62.96

Germany

wissenschaftlicherdienst@abcam.com | 030-896-779-154

Spain

soportecientifico@abcam.com | 91-114-65-60

Switzerland

technical@abcam.com

Deutsch: 043-501-64-24 | Français: 061-500-05-30

UK, EU and ROW

technical@abcam.com | +44(0)1223-696000

Canada

ca.technical@abcam.com | 877-749-8807

US and Latin America

us.technical@abcam.com | 888-772-2226

Asia Pacific

hk.technical@abcam.com | (852) 2603-6823

China

cn.technical@abcam.com | 400 921 0189 | +86 21 2070 0500

Japan

technical@abcam.co.jp | +81-(0)3-6231-0940

Singapore

sg.technical@abcam.com | 800 188-5244

Australia

au.technical@abcam.com | +61-(0)3-8652-1450

New Zealand

nz.technical@abc.com | +64-(0)9-909-7829