

Version 6 Last updated 19 December 2018

# ab155898 6-Phosphofructokinase Activity Assay Kit (Colorimetric)

For the rapid, sensitive and accurate measurement of 6-phosphofructokinase (PFK) activity in various samples.

[View kit datasheet: www.abcam.com/ab155898](http://www.abcam.com/ab155898)  
(use [www.abcam.cn/ab155898](http://www.abcam.cn/ab155898) for China, or [www.abcam.co.jp/ab155898](http://www.abcam.co.jp/ab155898) for Japan)

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

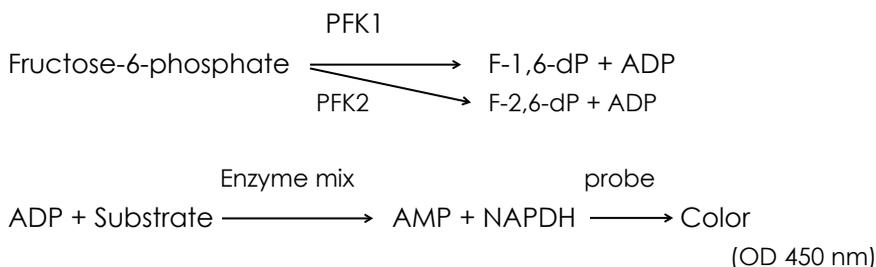
## Table of Contents

1. Overview	1
2. Protocol Summary	2
3. Precautions	3
4. Storage and Stability	3
5. Limitations	4
6. Materials Supplied	4
7. Materials Required, Not Supplied	5
8. Technical Hints	6
9. Reagent Preparation	7
10. Standard Preparation	9
11. Sample Preparation	10
12. Assay Procedure	12
13. Calculations	14
14. Typical Data	16
15. Quick Assay Procedure	18
16. Troubleshooting	19
17. Interferences	21
18. Notes	22

# 1. Overview

6-Phosphofructokinase Activity Assay Kit (Colorimetric) (ab155898) provides a simple and sensitive method to measure total 6-phosphofructokinase (PFK) activity in cell and tissue extracts. The assay is based in the conversion by PFK of fructose-6-phosphate to fructose-diphosphate, in presence of ATP. In the presence of a specific substrate and enzyme mix, The ADP produced from the reaction is converted to AMP and NADH, which in turn converts the colorless probe to a colored product with strong absorbance at OD 450 nm. The color intensity is proportional to the PFK activity present in the sample.

The assay can detect phosphofructokinase activity from as low as 1 mU PFK in solution.



Phosphofructokinase (PFK, EC 2.7.1.11) is a key glycolytic enzyme and plays a major regulatory role during glycolysis. This enzyme is present in bacteria, plants and animals. There are 2 types of phosphofructokinases: PFK1 and PFK2. In the presence of ATP, PFK1 catalyze the conversion of fructose-6-phosphate to fructose-1,6-diphosphate, whereas PFK2 catalyzes the conversion to fructose-2,6-diphosphate. PFK1 has three major isoforms in mammals: PFK1-M (muscle), PFK1-L (liver) and PFK1-P (platelet). In humans, PFK deficiency causes glycogen storage disease, also called Tarui's disease, which is characterized by exercise-induced muscle weakness and cramps. On the other hand, increased PFK activity contributes to cancer cell proliferation and tumorigenicity. Early detection of abnormal phosphofructokinase activity is crucial for diagnosis, prediction and therapeutic strategy.

## 2. Protocol Summary

Standard curve preparation



Sample preparation



Add reaction mix



Measure absorbance (OD450 nm) for 20-60 minutes at 37°C  
in kinetic mode\*

*\*For kinetic mode detection, incubation time given in this summary is for guidance only*

### 3. Precautions

**Please read these instructions carefully prior to beginning the assay.**

- All kit components have been formulated and quality control tested to function successfully as a kit.
- We understand that, occasionally, experimental protocols might need to be modified to meet unique experimental circumstances. However, we cannot guarantee the performance of the product outside the conditions detailed in this protocol booklet.
- Reagents should be treated as possible mutagens and should be handled with care and disposed of properly. Please review the Safety Datasheet (SDS) provided with the product for information on the specific components.
- Observe good laboratory practices. Gloves, lab coat, and protective eyewear should always be worn. Never pipet by mouth. Do not eat, drink or smoke in the laboratory areas.
- All biological materials should be treated as potentially hazardous and handled as such. They should be disposed of in accordance with established safety procedures.

### 4. Storage and Stability

**Store kit at -20°C in the dark immediately upon receipt. Kit has a storage time of 1 year from receipt, providing components have not been reconstituted.**

Refer to list of materials supplied for storage conditions of individual components. Observe the storage conditions for individual prepared components in the Materials Supplied section.

Aliquot components in working volumes before storing at the recommended temperature.

**Δ Note:** Reconstituted components are stable for 2 months.

## 5. Limitations

- Assay kit intended for research use only. Not for use in diagnostic procedures.
- Do not mix or substitute reagents or materials from other kit lots or vendors. Kits are QC tested as a set of components and performance cannot be guaranteed if utilized separately or substituted.

## 6. Materials Supplied

Item	Quantity	Storage temperature (before prep)	Storage temperature (after prep)
Assay Buffer	27 mL	-20°C	-20°C
PFK Substrate (lyophilized)	1 vial	-20°C	-20°C
ATP (lyophilized)	1 vial	-20°C	-20°C
PFK Enzyme Mix (lyophilized)	1 vial	-20°C	-20°C
PFK Developer (lyophilized)	1 vial	-20°C	-20°C
NADH Standard (lyophilized)	1 vial	-20°C	-20°C
Positive Control (lyophilized)	1 vial	-20°C	-20°C

## 7. Materials Required, Not Supplied

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

- Microplate reader capable of measuring absorbance at OD 450 nm
- Double distilled water (ddH<sub>2</sub>O)
- PBS
- Pipettes and pipette tips, including multi-channel pipette
- Assorted glassware for the preparation of reagents and buffer solutions
- Tubes for the preparation of reagents and buffer solutions
- 96-well plate with clear flat bottom
- Dounce homogenizer (if using tissue)
- (Optional) BCA protein assay kit (reducing agent compatible): we recommend using BCA protein assay kit reducing agent compatible (microplate) (ab207003)
- (Optional) 10 kD Spin Column (ab93349) – to filter sample lysates if background noise is very high

## 8. Technical Hints

- **This kit is sold based on number of tests. A “test” simply refers to a single assay well. The number of wells that contain sample, control or standard will vary by product. Review the protocol completely to confirm this kit meets your requirements. Please contact our Technical Support staff with any questions.**
- Selected components in this kit are supplied in surplus amount to account for additional dilutions, evaporation, or instrumentation settings where higher volumes are required. They should be disposed of in accordance with established safety procedures.
- Avoid foaming or bubbles when mixing or reconstituting components.
- Avoid cross contamination of samples or reagents by changing tips between sample, standard and reagent additions.
- Ensure plates are properly sealed or covered during incubation steps.
- Ensure all reagents and solutions are at the appropriate temperature before starting the assay.
- Samples generating values that are greater than the most concentrated standard should be further diluted in the appropriate sample dilution buffer.
- Make sure all necessary equipment is switched on and set at the appropriate temperature.

## 9. Reagent Preparation

Briefly centrifuge small vials at low speed prior to opening.

### 9.1 PFK Assay Buffer (50 mL):

Ready to use as supplied. Equilibrate to room temperature before use. Store at -20°C.

### 9.2 PFK Substrate (lyophilized, 1 vial):

Reconstitute the PFK substrate in 220 µL Assay Buffer. Aliquot so that you have enough to perform the desired number of assays. Store at -20°C. Avoid freeze/thaw cycles. Once reconstituted, use within two months. Keep on ice while in use.

### 9.3 ATP (lyophilized, 1 vial):

Reconstitute ATP in 220 µL ddH<sub>2</sub>O. Aliquot so that you have enough to perform the desired number of assays. Store at -20°C. Avoid freeze/thaw cycles. Once reconstituted, use within two months. Keep on ice while in use.

### 9.4 PFK Enzyme Mix (lyophilized, 1 vial):

Reconstitute PFK enzyme mix in 220 µL Assay Buffer. Aliquot so that you have enough to perform the desired number of assays. Store at -20°C. Avoid freeze/thaw cycles. Once reconstituted, use within two months. Keep on ice while in use.

### 9.5 PFK Developer (lyophilized, 1 vial):

Reconstitute PFK developer in 220 µL ddH<sub>2</sub>O. Aliquot so that you have enough to perform the desired number of assays. Store at -20°C. Avoid freeze/thaw cycles. Once reconstituted, use within two months. Keep on ice while in use.

### 9.6 NADPH Standard (lyophilized, 1 vial):

Reconstitute NADPH Standard in 40 µL Assay Buffer to generate 10 mM NADPH Standard solution. Aliquot standard so that you have enough volume to perform the desired number of assays. Store at -20°C. Use within 2 months. Keep on ice while in use.

### 9.7 Positive control (lyophilized, 1 vial):

Reconstitute positive control in 100  $\mu$ L Assay Buffer. Aliquot so that you have enough to perform the desired number of assays. Store at  $-20^{\circ}\text{C}$ . Avoid freeze/thaw cycles. Once reconstituted, use within two months. Keep on ice while in use.

## 10. Standard Preparation

- Always prepare a fresh set of standards for every use.
- Discard working standard dilutions after use as they do not store well.

**10.1** Prepare a 1 mM NADPH working standard solution (1:10 dilution) by adding 10  $\mu\text{L}$  10 mM NADPH Standard to 90  $\mu\text{L}$  Assay Buffer.

**10.2** Using 0.2 mM NADPH working standard, prepare standard curve dilution as described in the table in a microplate or microcentrifuge tubes:

Standard #	NADPH 1 mM Standard ( $\mu\text{L}$ )	Assay Buffer ( $\mu\text{L}$ )	Final volume standard in well ( $\mu\text{L}$ )	End amount NADPH in well (nmol/well)
1	0	150	50	0
2	6	144	50	2
3	12	138	50	4
4	18	132	50	6
5	24	126	50	8
6	30	120	50	10

Each dilution has enough amount of standard to set up duplicate readings (2 x 50  $\mu\text{L}$ ).

## 11. Sample Preparation

### General sample information:

- We recommend performing several dilutions of your sample to ensure the readings are within the standard value range.
- We recommend that you use fresh samples. If you cannot perform the assay at the same time, we suggest that you snap freeze your samples in liquid nitrogen upon extraction and store them immediately at  $-80^{\circ}\text{C}$ . When you are ready to test your samples, thaw them on ice and proceed with the Sample Preparation step. Be aware however that this might affect the stability of your samples and the readings can be lower than expected.

### 11.1 Cell lysates:

- 11.1.1 Harvest the amount of cells necessary for each assay (initial recommendation:  $2 \times 10^6$  cells).
- 11.1.2 Wash cells with cold PBS.
- 11.1.3 Resuspend cells in 100  $\mu\text{L}$  ice cold Assay buffer (containing protease inhibitors).
- 11.1.4 Homogenize cells quickly by pipetting up and down a few times.
- 11.1.5 Keep on ice for 10 minutes.
- 11.1.6 Centrifuge sample for 5 minutes at  $4^{\circ}\text{C}$  at 12,000  $\times g$  using a cold microcentrifuge to remove any insoluble material.
- 11.1.7 Collect supernatant and transfer to a new tube.
- 11.1.8 Keep on ice.
- 11.1.9 Optional: measure protein amount in the sample. Initial recommendation for reaction: 100  $\mu\text{g}$  protein/well.

### 11.2 Tissue lysates:

- 11.2.1 Harvest the amount of tissue necessary for each assay (initial recommendation  $\sim 20$  mg).
- 11.2.2 Wash tissue with cold PBS.
- 11.2.3 Homogenize tissue in 100  $\mu\text{L}$  of ice cold Assay Buffer (containing protease inhibitors) with a Dounce homogenizer sitting on ice, with 10 – 15 passes.
- 11.2.4 Centrifuge sample for 5 minutes at  $4^{\circ}\text{C}$  at 12,000  $\times g$  using a cold microcentrifuge to remove any insoluble material.
- 11.2.5 Collect supernatant and transfer to a new tube.

11.2.6 Keep on ice.

11.2.7 Optional: measure protein amount in the sample. Initial recommendation for reaction: 100 µg protein/well.

**Δ Note:** if sample background control wells show very high signal, you might need to remove interfering molecules with a 10 kD Spin Column (ab93349).

- Prewet the spin column with ddH<sub>2</sub>O and spin down ddH<sub>2</sub>O for 2 minutes at 10,000 x *g* at 4°C in a cold microcentrifuge. Remove ddH<sub>2</sub>O from upper and bottom reservoirs.
- Add 100 µL sample lysate and spin down for 10 minutes at 10,000 x *g* at 4°C. Discard filtrate and collect upper fraction.

**Δ Note:** We suggest using different volumes of sample to ensure readings are within the standard curve range.

## 12. Assay Procedure

- Equilibrate all materials and prepared reagents to room temperature prior to use.
- We recommend that you assay all standards, controls and samples in duplicate.
- Prepare all reagents, working standards, and samples as directed in the previous sections.
- A positive control is provided to ensure the assay is working correctly. It shouldn't be used as standard or to extrapolate enzyme activity from the sample.
- If sample background control wells show high signal, we recommend performing an additional filtration step with our 10 kD Spin Column (ab93349) to remove interfering molecules (See Section 11 for more details).

**Δ Note:** Small molecules such as glucose, ADP or NADPH present in cell or tissue extracts can generate background in this assay. We recommend that you set up Sample Background Controls to control for background noise.

### 12.1 Reaction wells set up:

- Standard wells = 50  $\mu$ L standard dilutions.
- Sample wells = 1-50  $\mu$ L samples (adjust volume to 50  $\mu$ L/well with Assay Buffer).

**Δ Note:** we recommend testing 100  $\mu$ g protein per well.

- Sample Background Control wells = 1-50  $\mu$ L samples (adjust volume to 50  $\mu$ L/well with Assay Buffer).

**Δ Note:** we recommend testing 100  $\mu$ g protein per well.

- Positive control wells = 10-20  $\mu$ L PFK positive control (adjust volume to 50  $\mu$ L/well with Assay Buffer).

## 12.2 PFK Reaction mix:

12.2.1 Prepare 50  $\mu\text{L}$  of PFK Reaction Mix and Background Mix for each reaction. Prepare a master mix to ensure consistency.

Component	Reaction Mix ( $\mu\text{L}$ )	Background Reaction Mix ( $\mu\text{L}$ )
Assay Buffer	42	44
PFK Enzyme Mix	2	2
PFK Developer	2	2
ATP	2	2
PFK Substrate	2	0

12.2.2 Add 50  $\mu\text{L}$  of Reaction Mix into each standard, positive control and sample wells.

12.2.3 Add 50  $\mu\text{L}$  of Background Reaction Mix into the background control sample wells.

12.2.4 Mix thoroughly.

### 12.3 Measurement:

12.3.1 Measure output at OD 450 nm on a microplate reader in kinetic mode for at least 20-60 minutes at 37°C protected from light.

**$\Delta$  Note:** Incubation time depends on the PFK activity in the samples. We recommend measuring OD in a kinetic mode, and choosing two time points (T1 and T2) to calculate the PFK activity of the samples. For standard curve, do not subtract A2 from A1. Standard curve can also be read in end point mode (i.e. at the end of incubation time).

## 13. Calculations

- Samples producing signals greater than that of the highest standard should be further diluted in appropriate buffer and reanalyzed, then multiply the concentration found by the appropriate dilution factor.
- Use only the linear rate for calculation.

### 13.1 Standard curve calculation:

- 13.1.1 Subtract the mean absorbance value of the blank (Standard #1) from all standard and sample readings. This is the corrected absorbance.
- 13.1.2 Average the duplicate reading for each standard.
- 13.1.3 Plot standard curve readings and draw the line of the best fit to construct the standard curve (most plate reader software or Excel can do this step). Calculate the trend line equation based on your standard curve data (use the equation that provides the most accurate fit).

### 13.2 Measurement of PFK activity in the sample:

- 13.2.1 For all reaction wells (including background control samples), choose two time points (T1 and T2) in the linear phase of the reaction progress curves and obtain the corresponding OD values at those points (OD1 and OD2)
- 13.2.2 Calculate  $\Delta OD$  for sample as follows:

$$\Delta OD_{450nm} = A2 - A1$$

- 13.2.3 Determine the background corrected change in fluorescence intensity for each well of sample by subtracting the  $\Delta OD$  value of the background control (BC).
- 13.2.4 PFK activity (nmol/min/mL or mU/mL) in the test samples is calculated as:

$$PFK \text{ Activity} = \left( \frac{B}{\Delta T \times V} \right) * D$$

Where:

B = amount of NADPH in sample well calculated from standard curve (nmol).

$\Delta T$  = linear phase reaction time T2 – T1 (minutes).

V = original sample volume added into the reaction well (mL).

D = sample dilution factor.

PFK activity can also be expressed as mU/ $\mu$ g of total protein in the sample.

Unit definition:

1 Unit PFK activity = amount of PFK that will generate 1.0  $\mu$ mol of NADPH per minute at pH 7.4 at 37°C.

## 14. Typical Data

Typical standard curve – data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.

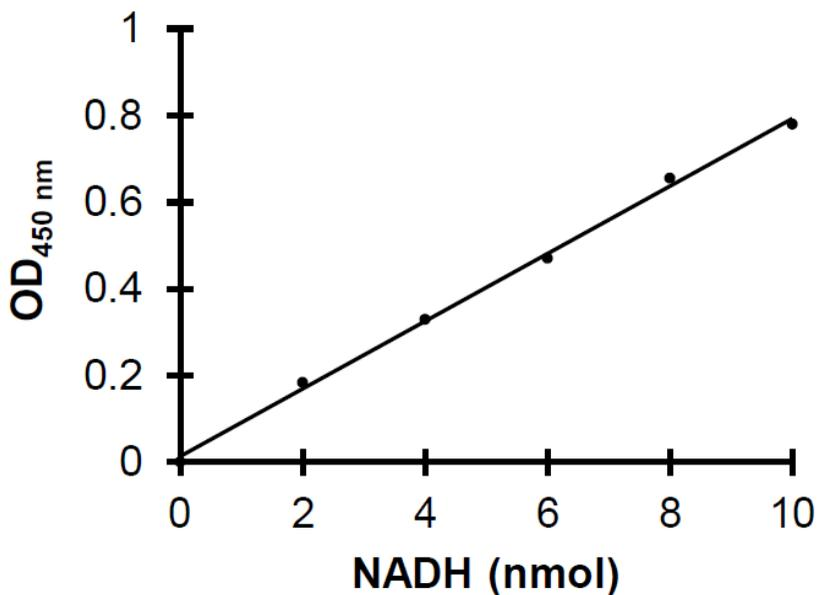
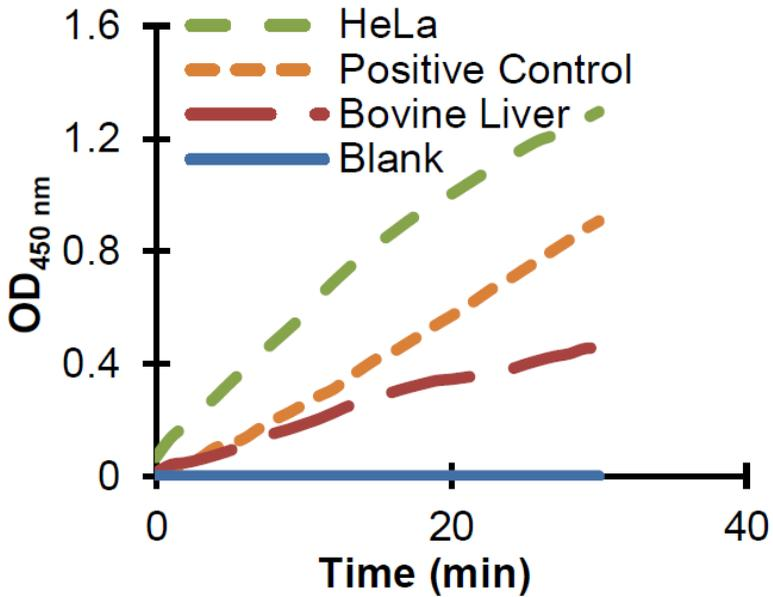


Figure 1. Typical NADH standard calibration curve.



**Figure 2.** Kinetic curves showing PFK activity detection in HeLa cell lysate (green), bovine liver lysate (red) and positive control included in the kit (orange). Blank (Standard #1) is shown in blue.

## 15. Quick Assay Procedure

**Δ Note:** this procedure is provided as a quick reference for experienced users. Follow the detailed procedure when performing the assay for the first time.

- Prepare reagents and aliquot; get equipment ready.
- Prepare PFK standard dilution [2 – 10 nmol/well].
- Prepare samples in optimal dilutions to fit standard curve readings.
- Set up plate in duplicate for standard (50  $\mu$ L), positive control (50  $\mu$ L), samples (50  $\mu$ L) and background sample control wells (50  $\mu$ L).
- Prepare a master mix for PFK Reaction Mix and (if appropriate) a master mix for Background Reaction Mix:

Component	Reaction Mix ( $\mu$ L)	Background Reaction Mix ( $\mu$ L)
Assay Buffer	42	44
PFK Enzyme Mix	2	2
PFK Developer	2	2
ATP	2	2
PFK Substrate	2	0

- Add 50  $\mu$ L Reaction to standard, positive control and sample wells.
- Add 50  $\mu$ L Background Reaction Mix to Sample Background control wells.
- Measure absorbance immediately at OD = 450 nm in a kinetic mode for 20-60 minutes at 37 °C protected from light.

## 16. Troubleshooting

Problem	Reason	Solution
<b>Assay not working</b>	Use of ice-cold buffer	Buffers must be at assay temperature
	Plate read at incorrect wavelength	Check the wavelength and filter settings of instrument
	Use of a different microplate	Colorimetric: clear plates Fluorometric: black wells/clear bottom plates Luminometric: white wells/clear bottom plates
<b>Sample with erratic readings</b>	Cells/tissue samples not homogenized completely	Use Dounce homogenizer, increase number of strokes
	Samples used after multiple free/ thaw cycles	Aliquot and freeze samples if needed to use multiple times
	Use of old or inappropriately stored samples	Use fresh samples or store at - 80°C (after snap freeze in liquid nitrogen) till use
	Presence of interfering substance in the sample	Check protocol for interfering substances; deproteinize samples
<b>Lower/higher readings in samples and standards</b>	Improperly thawed components	Thaw all components completely and mix gently before use
	Allowing reagents to sit for extended times on ice	Always thaw and prepare fresh reaction mix before use
	Incorrect incubation times or temperatures	Verify correct incubation times and temperatures in protocol

<b>Problem</b>	<b>Reason</b>	<b>Solution</b>
<b>Standard readings do not follow a linear pattern</b>	Pipetting errors in standard or reaction mix	Avoid pipetting small volumes (< 5 $\mu$ L) and prepare a master mix whenever possible
	Air bubbles formed in well	Pipette gently against the wall of the tubes
	Standard stock is at incorrect concentration	Always refer to dilutions described in the protocol
<b>Unanticipated results</b>	Measured at incorrect wavelength	Check equipment and filter setting
	Samples contain interfering substances	Troubleshoot if it interferes with the kit
	Sample readings above/ below the linear range	Concentrate/ Dilute sample so it is within the linear range

## 17. Interferences

These chemical or biological materials will cause interferences in this assay causing compromised results or complete failure:

- Small molecules such as glucose, NADH or ADP present in the sample: we recommend setting up a background control well to correct for noise interference.
- RIPA buffer: SDS may interfere with the enzymatic reaction.

## 18. Notes







## Technical Support

Copyright © 2017 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@abcam.com | +64-(0)9-909-7829