

Version 2 Last updated 28 April 2021

ab234624 Collagenase (Collagen Degradation/Zymograph y) Assay Kit (Fluorometric)

For the measurement of collagenase activity in biological samples.

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	3
5. Reagent Preparation	4
6. Standard Preparation	5
7. Sample Preparation	6
8. Assay Procedure	8
9. Data Analysis	9
10. Typical Data	10
11. Notes	12

1. Overview

Collagenase (Collagen Degradation/Zymography) Assay Kit (Fluorometric) (ab234624) employs a highly quenched collagen substrate which upon cleavage by a suitable collagenase releases a fluorophore, which can be easily quantified using a conventional microplate reader. This method is substrate-specific, simple, fast, high-throughput adaptable and amenable to the sensitive detection of collagenase activity (as low as 0.6 mCDU for bacterial collagenase) in biological samples.

Prepare samples



Prepare FITC Standards



Prepare Collagenase Substrate Mix



Add Substrate Mix, Samples and Positive Controls into wells and measure fluorescence in kinetic mode Ex/Em 490/520 nM

2. Materials Supplied and Storage

Store kit at -20°C in the dark immediately on receipt and check below for storage for individual components. Kit can be stored for 1 year from receipt, if components have not been reconstituted.

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

Avoid repeated freeze-thaws of reagents.

Item	Quantity	Storage temperature (before prep)	Storage temperature (after prep)
Cell Lysis Buffer	25 mL	-20°C	-20°C
Collagenase Assay Buffer	25 mL	-20°C	-20°C
Collagenase Substrate	1 vial	-20°C	-20°C
Enzyme Positive Control	100 µL	-20°C	-20°C
FITC Standard (5 mM)	10 µL	-20°C	-20°C

3. 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 fluorescence at Ex/Em = 490/520 nm
- 96 well plate with clear flat bottom, preferably black (for fluorometric assay)
- Dounce homogenizer

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. Reagent Preparation

Briefly centrifuge small vials at low speed prior to opening.

5.1 Collagenase Assay Buffer

Ready to use as supplied. Bring to room temperature before use.

5.2 Cell Lysis Buffer

Ready to use as supplied.

5.3 Enzyme Positive Control

Ready to use as supplied. Thaw on ice before use.

5.4 Collagenase Substrate

1. Reconstitute in 220 μL of deionized water.
2. Mix well by pipetting up and down. Vortex if necessary.
3. Unused substrate can be stored at -20°C by covering it with aluminum foil or transferring it to an amber vial.

5.5 FITC Standard (5 mM)

Ready to use as supplied.

6. Standard Preparation

- Always prepare a fresh set of standards for every use.
 - Discard working standard dilutions after use as they do not store well.
1. Prepare a 50 μM FITC standard by diluting 2 μL of 5 mM FITC Standard with 198 μL of Collagenase Assay Buffer. Mix well by pipetting up and down, vortex vigorously for 30 seconds.
 2. Using the 50 μM standard, prepare standard curve dilution as described in the table in a microplate or microcentrifuge tubes:

Standard #	FITC Standard (μL)	Assay Buffer (μL)	Final volume standard in well (μL)	End amount of FITC standard in well (pmol/well)
1	0	200	100	0
2	4	196	100	100
3	8	192	100	200
4	12	188	100	300
5	16	184	100	400
6	20	180	100	500

Each dilution has enough standard to set up duplicate readings (2 x 100 μL).

3. Mix well and measure the fluorescence at Ex/Em 490/520 nm in end-point mode at 37 $^{\circ}\text{C}$.

7. 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 for the most reproducible assay. 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°C . When you are ready to test your samples, thaw them on ice. Be aware, however, that this might affect the stability of your samples and the readings can be lower than expected. Avoid multiple freeze-thaws.

7.1 Fresh/frozen tissue and cells:

1. Homogenize fresh or frozen tissue (~5-10 mg) or cells ($1-2 \times 10^6$) with 100 μL Cell Lysis Buffer and incubate on ice for 5 minutes.
2. Centrifuge the homogenate at $16,000 \times g$, 4°C for 10 minutes.
3. Transfer the clarified supernatant to a fresh pre-chilled tube and keep on ice.
4. Measure the amount of protein in the lysate or purified enzyme using BCA protein assay kit reducing agent compatible (microplate) (ab207003).
5. Add 1-50 μL of lysate or purified enzyme into desired well(s) in a white 96-well plate. If necessary, dilute the lysate with Collagenase Assay buffer.
6. For the Positive Control, dilute 2 μL of Enzyme Positive Control with 18 μL of Collagenase Assay Buffer and use 1-10 μL /well.
7. Adjust the volume of Samples and Positive Control to 50 μL /well with Collagenase Assay Buffer.

Δ Note: *The kit is designed to work with active Collagenase enzymes only. If the sample contains inactive zymogen forms of collagenase, it can be activated with p-aminophenylmercuric acetate (APMA) or other activators. The conditions for activation of each enzyme should be determined empirically by following an appropriate testing protocol (Shapiro et. al., J. Bio. Chem. 1995, 270 (11), 6351-6356).*

Δ Note: *To induce higher collagenase expression, cells can be grown in the presence of Phorbol myristate acetate (10-50 ng/ml), lysed and tested directly in the assay (Hersh et. al., Biochem., 1986, 25 (17), 4750-4757).*

Δ Note: For unknown samples, we suggest doing pilot experiment and testing 3-5 different amounts of samples to ensure the readings are within the Standard Curve range.

8. Assay Procedure

- Equilibrate all materials and prepared reagents to room temperature just prior to use and gently agitate, apart from Positive Control, keep this on ice.
- Assay all standards, controls and samples in duplicate.

Δ Note: *If you suspect your samples contain substance that can generate background, set up Sample Background Controls to correct for background noise.*

8.1 Collagenase Substrate Mix:

1. Prepare 50 μL of Collagenase Substrate Mix per well. Prepare a master mix to ensure consistency.
2. Dissolve the Substrate Mix by vigorous vortexing.

Component	Substrate Mix (μL)
Collagenase Assay Buffer	48
Collagenase Substrate	2

3. Add 50 μL of Substrate Mix into each Sample and Positive Control well.
4. Mix well and measure the fluorescence at Ex/Em 490/520 nm in kinetic mode at 37 °C for 1-2 hours.

Δ Note: *Do not add Substrate Mix to the sample Background Control and Standard wells.*

9. Data Analysis

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.

1. Choose two time points (t1 and t2) where the corresponding RFUs (RFU1 and RFU2) are in a linear range.
2. Calculate ΔRFU and Δt and obtain $\Delta\text{RFU}/\Delta t$ as RFU/minute for each Sample including Background Control.
3. Subtract the value of RFU/minute of Background Control from each Sample to obtain net RFU/minute (B).
4. Obtain change in the RFU (ΔRFU) by subtracting fluorescence of the 0 Standard Controls from those containing all standards.
5. Plot the ΔRFU against pmol of FITC Standard.
6. Apply the corrected sample RFU reading to the standard curve to get Collagenase (B) amount in the sample wells.
7. The plot should be linear; determine the slope A ($\Delta\text{RFU}/\text{pmol}$) of the curve.
8. Using RFU/min of each Sample, calculate Sample Collagenase activity using following equation:

$$\text{Collagenase activity } \left(X, \frac{\text{U}}{\text{ml}} \right) = \frac{\text{B} \times 1000}{\text{A} \times \text{C}} * \text{D}$$

$$\text{Collagenase activity } \left(\frac{\text{U}}{\text{mg}} \right) = X/P$$

Where:

B = Sample Collagenase Activity as calculated (RFU/minute)

A = Slope of the FITC standard curve ($\Delta\text{RFU}/\text{pmol}$)

C = μl of Sample used in the assay

D = sample dilution factor if sample is diluted to fit within the standard curve range (prior to reaction well set up).

P = Protein concentration in the lysate (mg/mL)

1000 = Conversion Factor ($1000 \mu\text{L} \equiv 1 \text{ mL}$)

Unit Definition: 1 U is the amount of Collagenase required to cleave the Collagenase Substrate and release 1 pmol of Fluorescein per minute under the conditions of the assay.

10. Typical Data

Data provided for demonstration purposes only.

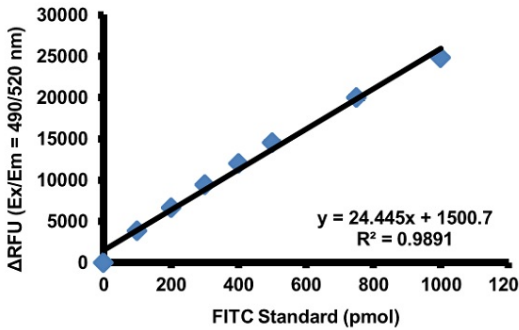


Figure 1. Example Standard Curve.

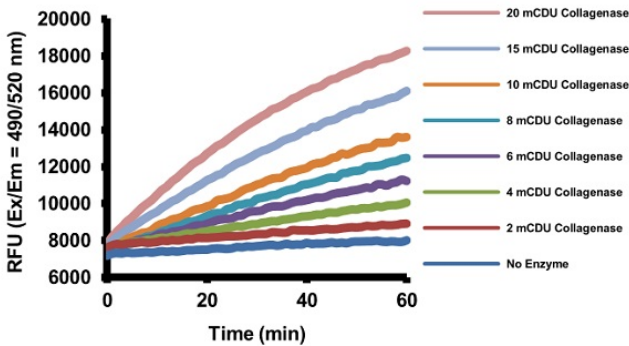


Figure 2. Collagenase activity with different amounts of Enzyme Positive Control.

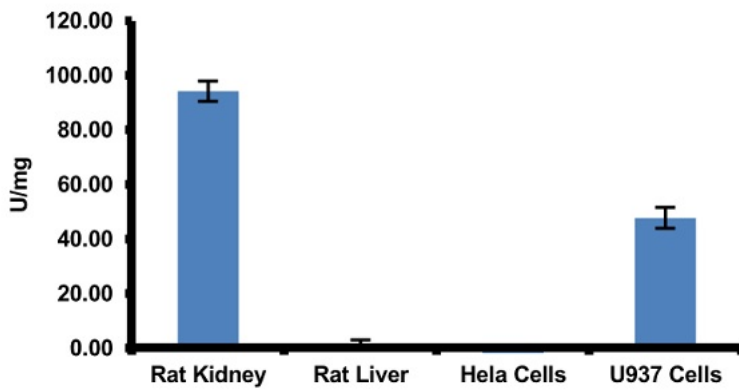


Figure 3. Collagenase activity in rat kidney, liver lysates along with HeLa and U937 cell lysates. n=3.

11. Notes

Technical Support

Copyright © 2021 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 | +86-21-5110-5938 | 400-628-6880

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