

## Product datasheet

# Gelatin Degradation Assay Kit (Alternative to Zymography) ab234057

[1 References](#) [3 Images](#)

### Overview

<b>Product name</b>	Gelatin Degradation Assay Kit (Alternative to Zymography)
<b>Detection method</b>	Fluorescent
<b>Sample type</b>	Tissue, Cell Lysate
<b>Assay type</b>	Quantitative
<b>Species reactivity</b>	<b>Reacts with:</b> Mammals, Other species
<b>Product overview</b>	Gelatin Degradation Assay Kit (Alternative to Zymography) (ab234057) utilizes a hybrid approach for the detection of gelatinase activity by employing a highly quenched gelatin substrate which upon cleavage by a suitable gelatinase 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 gelatinase activity (as low as 0.06 mCDU for bacterial collagenase) in biological samples.

<b>Notes</b>	<p>This product is manufactured by BioVision, an Abcam company and was previously called K444 Gelatinase (Gelatin Degradation/Zymography) Assay Kit (Fluorometric). K444-100 is the same size as the 100 test size of ab234057.</p> <p>Gelatinases are a type of matrix zinc-dependent metalloproteases (MMPs) that degrade gelatins and a variety of other extracellular matrix proteins. These enzymes are synthesized as latent zymogens that are activated by proteolysis and inhibited by tissue inhibitors of metalloproteases (TIMPs). Two mammalian gelatinases, Gelatinase A (MMP-2) and Gelatinase B (MMP-9), are critical for basement membrane degradation and are highly upregulated in variety of tumor cells. Gelatinase activity is usually detected by small peptide-based activity assays which may suffer from lack of substrate specificity. Other methods for gelatinase activity include gelatin Zymography where samples are electrophoresed on a gelatin-containing SDS-PAGE, and further renatured in a suitable buffer for 12-16 h. The zymogram is subsequently stained, and areas of digestion appear as clear bands against a darkly stained background where the substrate has been degraded by the enzyme. Such methods are laborious, time-consuming and may lead to the loss of enzymatic activity as renaturation may not be completely reversible.</p>
--------------	---

<b>Platform</b>	Microplate reader
-----------------	-------------------

### Properties

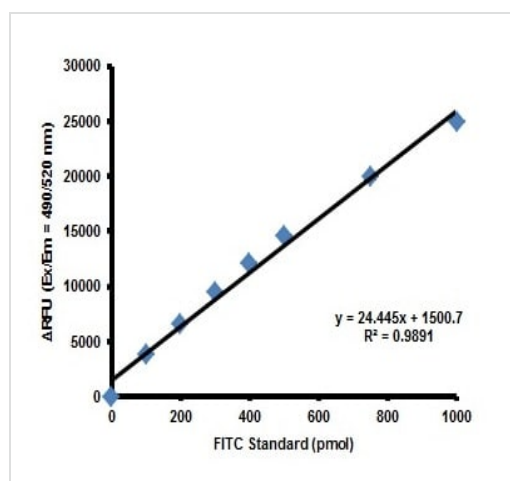
<b>Storage instructions</b>	Store at -20°C. Please refer to protocols.
-----------------------------	--

Components	100 tests
Cell Lysis Buffer	1 x 25ml
Enzyme Positive Control	1 x 100µl
FITC Standard (5 mM)	1 x 10µl
Gelatinase Assay Buffer	1 x 25ml
Gelatinase Substrate (FITC)	1 vial

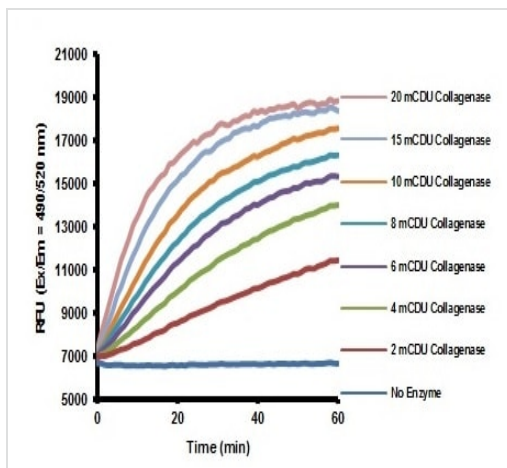
## Relevance

Gelatin is denatured collagen composed of heterogeneous mixture of water-soluble proteins of high average molecular weights. Below 35-40°C gelatin swells and absorbs 5-10 times its weight of water to form a gel. Gel strength and viscosity gradually weaken upon prolonged heating in solution above 40°C.

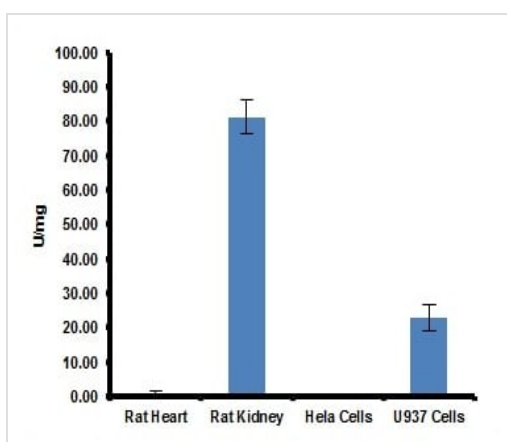
## Images



FITC Standard Curve.



Gelatinase activity with different amounts of Enzyme  
Positive Control.



Gelatinase activity in rat heart and kidney lysates  
along with HeLa and U937 cell lysates.

**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