

# Collagenase Activity Assay Kit (Colorimetric) ab196999

[3 References](#) [2 Images](#)

### Overview

<b>Product name</b>	Collagenase Activity Assay Kit (Colorimetric)
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Cell Lysate, Purified protein
<b>Assay type</b>	Enzyme activity
<b>Sensitivity</b>	> 20 $\mu$ U
<b>Assay time</b>	0h 25m
<b>Assay duration</b>	Multiple steps standard assay
<b>Species reactivity</b>	<b>Reacts with:</b> Bacteria
<b>Product overview</b>	Collagenase Activity Assay Kit (ab196999) provides a quick and easy way to determine activity of Collagenase. The assay measures collagenase activity using a synthetic peptide (FALGPA) that mimics the structure of collagen. It is suitable for measuring activity of bacterial collagenases such as from <i>Clostridium histolyticum</i> type I-XI. This assay can also be used for screening or characterization of collagenase inhibitors.

The limit of detection for this assay is 0.02 mU Collagenase.

Collagenase assay protocol summary:

- add samples and controls to wells
- add reaction mix
- analyze with microplate reader in kinetic mode for 5-15 min

**Notes** This product is manufactured by BioVision, an Abcam company and was previously called K792 Collagenase Activity Colorimetric Assay Kit. K792-100 is the same size as the 100 test size of ab196999.

Collagenase (EC 3.4.24.3) is an enzyme in the matrix metalloproteinase family that breaks down collagen, assisting in degradation of the extracellular matrix, which is a key step in the pathogenesis of bacteria. Collagen is an abundant structural protein present in the connective tissue of animals. Collagenase has been used clinically for the treatment of Dupuytren's contracture, an affliction characterized by a thickening of connective tissue.

**Platform** Microplate reader

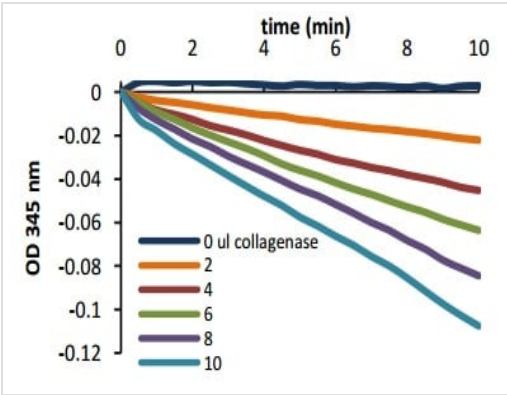
### Properties

Storage instructions

Store at -20°C. Please refer to protocols.

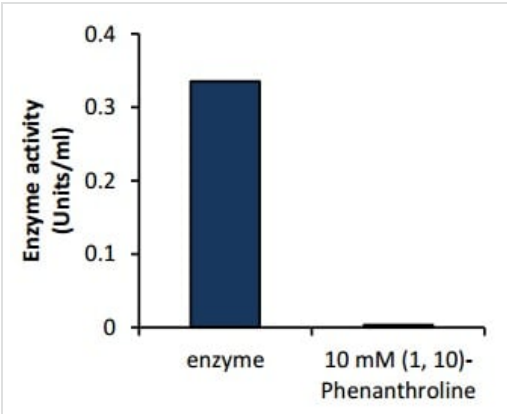
Components	100 tests
Collagenase	1 x 1ml
Collagenase Assay Buffer	1 x 20ml
Collagenase Substrate I	1 x 4ml
Inhibitor (1,10-Phenanthroline)	1 x 50µl

Images



Collagenase activity over 10 minutes.

Functional Studies - Collagenase Activity Assay Kit  
(Colorimetric) (ab196999)



Enzyme activity of collagenase and inhibition by 10 mM (1,10)-Phenanthroline. Assay was performed following the kit protocol.

Functional Studies - Collagenase Activity Assay Kit  
(Colorimetric) (ab196999)

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