

# Enhanced Apoptotic DNA Ladder Detection Kit ab66093

### Overview

|                         |  |
|-------------------------|--|
| <b>Product name</b>     | Enhanced Apoptotic DNA Ladder Detection Kit  |
| <b>Sample type</b>      | Tissue, Adherent cells, Suspension cells   |
| <b>Assay type</b>       | Direct   |
| <b>Assay time</b>       | 1h 30m   |
| <b>Product overview</b> | Enhanced Apoptotic DNA Ladder Detection Kit (ab66093) provides an easy and sensitive solution for detecting DNA fragmentation in apoptotic cells. It can isolate small fragmented DNA from cells in only 90 minutes, whereas other similar products in the market require 1-2 days. No DNA extraction or additional columns required. DNA fragments can be easily visualized by agarose gel electrophoresis stained with a highly sensitive dye. |

|              |  |
|--------------|--|
| <b>Notes</b> | <p>This product is manufactured by BioVision, an Abcam company and was previously called K130 Enhanced Apoptotic DNA Ladder Detection Kit. K130-50 is the same size as the 50 test size of ab66093.</p> <p>For more apoptosis assays, review the <a href="#">apoptosis assay and apoptosis marker guide</a>.</p> |
|--------------|--|

### Properties

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

| Components                | 50 tests   |
|---------------------------|------------|
| Ammonium Acetate Solution | 1 x 250µl  |
| DNA Suspension Buffer     | 1 x 0.25ml |
| Enzyme B Mix              | 1 vial     |
| Staining Dye              | 1 x 50µl   |
| TE Lysis Buffer           | 1 x 1.8ml  |

|                  |   |
|------------------|---|
| <b>Relevance</b> | Internucleosomal DNA fragmentation is a hallmark of apoptosis in mammalian cells. |
|------------------|---|

**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