

# Apoptosis DNA Ladder Assay Kit ab66090

★★★★★ [1 Abreviews](#) [9 References](#) [1 Image](#)

### Overview

<b>Product name</b>	Apoptosis DNA Ladder Assay Kit
<b>Sample type</b>	Adherent cells, Suspension cells
<b>Assay type</b>	Direct
<b>Assay time</b>	1h 30m
<b>Product overview</b>	Apoptosis DNA Ladder Assay Kit (ab66090) 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. No DNA extraction or additional columns required.

Apoptosis DNA ladder assay protocol summary:

- pellet cells by spinning
- wash cells with PBS and spinning
- lyse cells with TE lysis buffer and pipetting
- add enzyme A solution and incubate for 10 min
- add enzyme B solution and incubate for 30 min
- add ammonium acetate solution, isopropanol and freeze for 10 min
- spin to precipitate DNA
- wash pellet with 70% ethanol, air-dry and dissolve in DNA suspension buffer
- run on agarose gel and visualize with DNA staining dye

**Notes** This product is manufactured by BioVision, an Abcam company and was previously called K120 Quick Apoptotic DNA Ladder Detection Kit. K120-50 is the same size as the 50 test size of ab66090.

For more apoptosis assays, review the [apoptosis assay and apoptosis marker guide](#).

### Properties

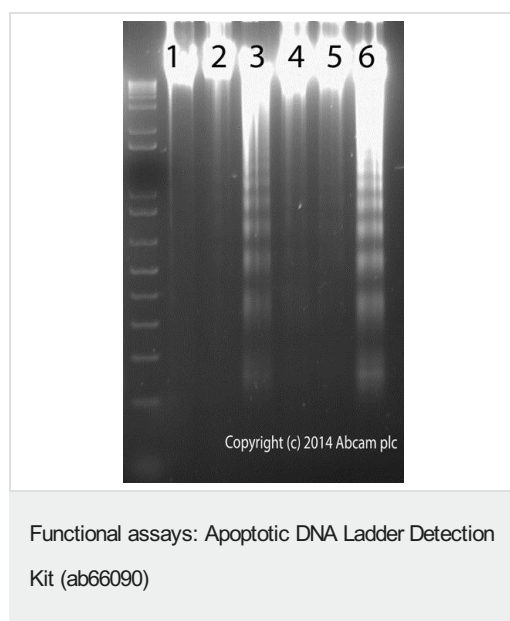
**Storage instructions** Store at -20°C. Please refer to protocols.

Components	50 tests
Ammonium Acetate/Glycogen	1 x 250µl
DNA Suspension Buffer	1 x 1.5ml

Components	50 tests
Enzyme A Solution	1 x 250µl
Enzyme B Mix	1 vial
Lysis Buffer III	1 x 1.8ml

**Relevance** Internucleosomal DNA fragmentation is a hallmark of apoptosis in mammalian cells.

## Images



Apoptosis was induced in 10e6 Jurkat cells by treatment with 10 µM Camptothecin (CPT) ([ab120115](#)) for 4 hours (2, 5) or 2 µM CPT for 24 hours (3, 6) whilst control cells were kept without treatment (1, 4). DNA was extracted and resuspended in 60 µL resuspension buffer. 10 µL (1-3) or 40 µL (4-6) were loaded onto the agarose gel.

**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