

Enhanced Apoptotic DNA Ladder Detection Kit ab66093

概述

产品名称	Enhanced Apoptotic DNA Ladder Detection试剂盒
样品类型	Tissue, Adherent cells, Suspension cells
检测类型	Direct
检测时间	1h 30m
产品概述	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.

说明	<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 apoptosis assay and apoptosis marker guide.</p>
----	--

性能

存放说明	Store at -20°C. Please refer to protocols.
------	--

组 件	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

相关性	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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors