

Product datasheet

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. |

Visit our [FAQs page](#) for tips and troubleshooting.

| | |
|-------|---|
| 说明 | Internucleosomal DNA fragmentation is a hallmark of apoptosis in mammalian cells. |
| 经测试应用 | 适用于: Functional Studies |

性能

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

| 组件 | 标识符 | 50 tests |
|---------------------------|--------|------------|
| Ammonium Acetate Solution | Yellow | 1 x 250µl |
| DNA Suspension Buffer | Green | 1 x 0.25ml |
| Enzyme B (Lyophilized) | Red | 1 vial |
| Staining Dye (10000X) | Blue | 1 x 50µl |
| TE Lysis Buffer | Purple | 1 x 1.8ml |

| | |
|-----|---|
| 相关性 | Internucleosomal DNA fragmentation is a hallmark of apoptosis in mammalian cells. |
|-----|---|

应用

Our [Abpromise guarantee](#) covers the use of **ab66093** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

| 应用 | Ab评论 | 说明 |
|--------------------|------|--|
| Functional Studies | | Use at an assay dependent concentration. |

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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 <http://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