

Apoptotic DNA Ladder Isolation Kit ab65627

6 References [画像数 2](#)

製品の概要

製品名	Apoptotic DNA Ladder Isolation Kit
サンプルの種類	Tissue, Adherent cells, Suspension cells
アッセイタイプ	Direct
製品の概要	Apoptotic DNA Ladder Isolation Kit (ab65627) provides an easy and sensitive solution for detecting DNA fragmentation in apoptotic cells and tissues. The new procedure selectively extracts DNA ladders without interference of intact genomic DNA in only 2-4 hours. As little as 5% of apoptotic cells in a sample can be visualized with this high sensitivity assay.

特記事項

This product is manufactured by BioVision, an Abcam company and was previously called K170 Apoptotic DNA Ladder Isolation Kit. K170-50 is the same size as the 50 test size of ab65627. For more apoptosis assays, review the [apoptosis assay and apoptosis marker guide](#).

製品の特性

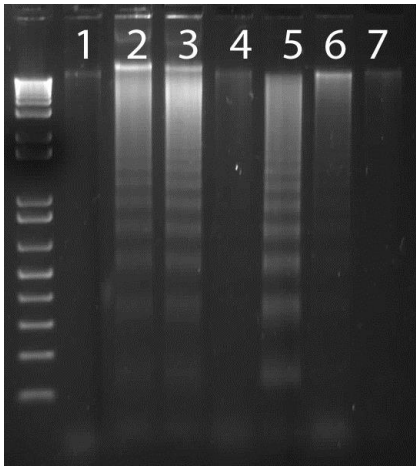
保存方法 Store at -20°C. Please refer to protocols.

内容	50 tests
Ammonium Acetate Solution	1 x 250µl
DNA Ladder Extraction Buffer	1 x 12.5ml
DNA Suspension Buffer	1 x 2ml
Enzyme A Solution	1 x 250µl
Enzyme B (Lyophilized)	1 vial

関連性

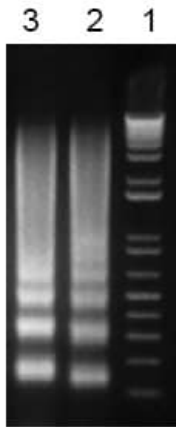
Internucleosomal DNA fragmentation is a hallmark of apoptosis in mammalian cells.

画像



Functional Assay: ab65627 Apoptotic DNA Ladder Isolation Kit

Apoptosis induced in Jurkat cells (1×10^6 cell \times mL⁻¹) by incubation for 24 hours with 1) Control medium; 2) 2 μ M Camptothecin (**ab120115**); 3) 10 μ M Camptothecin; 4) 10 μ M Camptothecin + 20 nM caspase inhibitor Z-VAD(OMe)-FMK (zVAD; **ab120487**); 5) 50 ng \times mL⁻¹ anti-Fas Ab (MBL), 6) 0.05 mM H₂O₂ (Sigma); 7) 1 mM H₂O₂. 5 μ L of DNA preparation loaded into each well.



Functional Studies - Apoptotic DNA Ladder Isolation Kit (ab65627)

DNA ladders were extracted from 2×10^6 Jurkat cells were treated with 2 μ M Camptothecin. The kit extracts laddered DNA only but not intact genomic DNA; therefore, more cells can be extracted and load on the agarose gel. Lane 1: 1 kb DNA ladder; Lane 2: 6 hour induction; Lane 3: 12 hour induction.

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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