

Product datasheet

Chromatin Extraction Kit ab117152

[38 References](#) [2 Images](#)

Overview

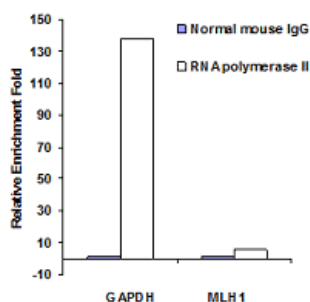
Product name	Chromatin Extraction Kit
Sample type	Tissue, Adherent cells, Suspension cells
Assay time	1h 00m
Product overview	Chromatin Extraction Kit (ab117152) is a complete set of optimized buffers and reagents for isolating chromatin or DNA-protein complex from mammalian cells or tissues in a simple and rapid format. Chromatin prepared by this kit can be used in a variety of chromatin immunoprecipitation methods. The isolated chromatin can also be used in other chromatin-related applications such as in vitro protein-DNA binding assays and nuclear enzyme assays.

Properties

Storage instructions Store at +4°C. Please refer to protocols.

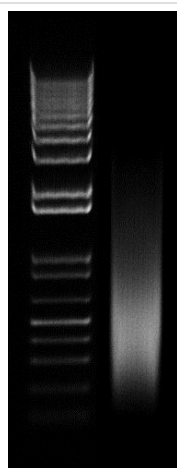
Components	100 tests
1000X Protease Inhibitor Cocktail	1 x 110µl
10X Lysis Buffer	1 x 11ml
Chromatin Buffer	1 x 11ml
Extraction Buffer	1 x 11ml

Images



ChIP analysis of RNA polymerase II enriched in GAPDH and MLH1 promoters with chromatin extract prepared from formaldehyde fixed colon cancer cells (2×10^5) using ab117152.

Functional Studies - Chromatin Extraction kit
(ab117152)



Chromatin was extracted from 3×10^6 mouse neuroblastoma cells, fixed for 10min at RT in 0.1% formaldehyde. Cells were sonicated before last centrifugation, and following centrifugation, cells were treated with proteinase K & RNase A. Chromatin yield was 4 μ g chromatin/ 10^6 cells. Lane 1: 1kb DNA ladder Lane 2: 300 - 650bp DNA fragments

Image courtesy of anonymous Abreview

Chromatin Extraction Kit (ab117152)

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