

Product datasheet

Nuclear Extraction Kit ab113474

★★★★★ 3 Abreviews 67 References 3 Images

Overview

Product name	Nuclear Extraction Kit
Sample type	Tissue, Adherent cells, Suspension cells
Assay time	1h 00m
Product overview	Nuclear Extraction Kit (ab113474) provides a simple and selective method along with all necessary reagents for nuclear protein extraction / nuclear protein fractionation in just 1 hour.

The extracts can then be used in western blotting, protein-DNA binding assays, nuclear enzyme assays or any other procedures requiring optimized nuclear proteins. The protocol is fast and easy-to-use, and isolates very abundant yields of nuclear extract from mammalian cells or tissue samples.

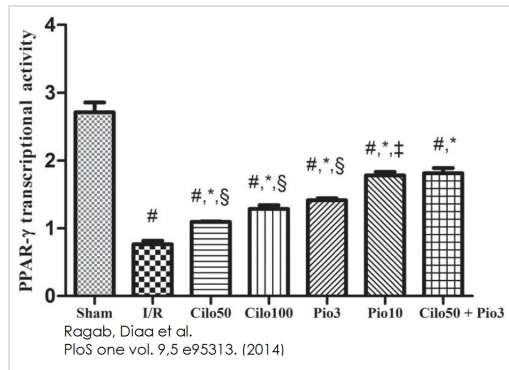
Not sure if this is the right product for you? Check out our [EpiSeeker Sample Preparation Guide](#) for help.

Compared to other kits that use conventional nuclear extraction / nuclear fractionation methods, the buffers included in ab113474 contain much lower amounts of salts (80% less than conventional kits) and no SDS, which allows much better retention of enzyme activity in the nuclear extracts.

Properties

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

Components	100 tests
1000X Protease Inhibitor Cocktail	1 x 110µl
10X Pre-Extraction Buffer	1 x 10ml
DTT Solution (1000X)	1 x 110µl
ENE2 (Extraction Buffer)	1 x 10ml



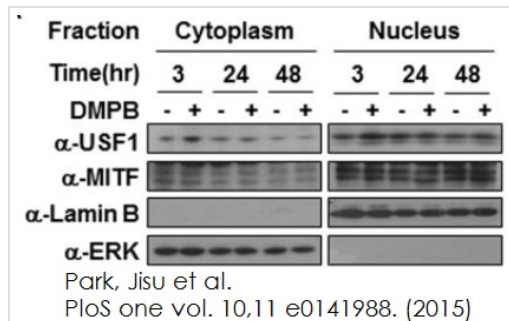
Functional Studies - Nuclear Extraction Kit

(ab113474)

Ragab, Diao et al., PloS one?vol. 9,5 e95313., Fig 6,
doi:10.1371/journal.pone.0095313

Effect of two drugs referred to as Cilo (50 and 100 mg/kg; Cilo50and Cilo100), and Pio (3 and 10 mg/kg; Pio3 and Pio10), and their combination (Cilo50and Pio3) on the PPAR-γ transcription activity in rats subjected to ischemia (45 min)/reperfusion (24 hrs).

Drugs were administered orally for 14 days then subjected to ischemia/reperfusion. Values are expressed as mean ± S.E.M (n = 6). Data are compared with sham operated control (#), I/R control (□), Cilo50 (○), Pio3 (△), and combination (§) pretreated groups (one-way ANOVA followed by Tukey Multiple Comparison Test) at P<0.05.

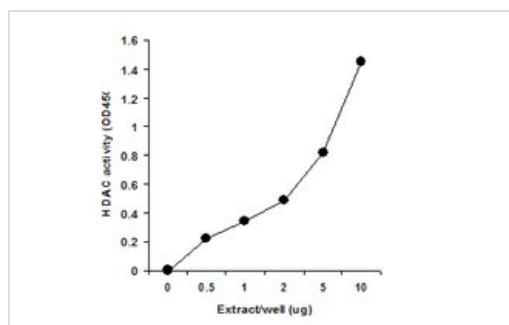


Functional Studies - Nuclear Extraction Kit

(ab113474)

Park, Jisu et al., PloS one?vol. 10,11 e0141988., Fig 4,
doi:10.1371/journal.pone.0141988

B16F10 cells were treated with 30 μM of DMPB for the indicated time periods. Cytoplasmic and nuclear fractions were isolated and analyzed by Western blotting.



Functional Studies - Nuclear Extraction Kit

(ab113474)

Nuclear extracts were prepared from MCF-7 cells and the activity of HDACs were measured using different amounts of the extract. The result shown in the figure demonstrates the ab113474's high specificity.

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