

## Product datasheet

# Nuclear Extraction Kit ab221978

### Overview

<b>Product name</b>	Nuclear Extraction Kit
<b>Sample type</b>	Tissue, Adherent cells, Suspension cells
<b>Product overview</b>	This Nuclear Extraction Kit is formulated for the quick and simple isolation of nuclear and cytoplasmic fractions from cultured cells and tissue homogenates that can be used successfully in our transcription factor assay kits.

### Notes

### Properties

**Storage instructions** Please refer to protocols.

Components	100 tests
Nonidet P-40 Assay Reagent (10%)	1 x 6ml
Nuclear Extraction Buffer (2X)	1 x 3ml
Nuclear Extraction Dithiothreitol (1M)	1 x 60µl
Nuclear Extraction Hypotonic Buffer (10X)	1 x 6ml
Nuclear Extraction PBS (10X)	1 x 100ml
Nuclear Extraction Phosphatase Inhibitors (50X)	2 x 5ml
Nuclear Extraction Protease Inhibitor Cocktail (100X)	1 x 650µl

**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