

Nuclear (Nucleic Acid-Free) Extraction Kit ab113477

3 References

概述

产品名称	核(Nucleic Acid-Free) Extraction试剂盒
样品类型	Tissue, Adherent cells, Suspension cells
检测时间	1h 00m
种属反应性	与反应: Mammals
产品概述	<p>Nuclear (Nucleic Acid-Free) Extraction Kit (ab113477) provides a simple and selective method for extracting nuclear proteins (nucleic-acid free) in just 1 hour for a variety of applications such as western blotting, protein-DNA binding assays, or any other procedures requiring optimized nucleic acid-free nuclear proteins.</p> <p>ab113477 can be used to extract nuclear proteins from mammalian cells and tissue samples free of nucleic acids attached.</p> <p>Not sure if this is the right product for you? Check out our Methods and tools to study histone modifications for help.</p>

性能

存放说明	Please refer to protocols.
------	----------------------------

组件	100 tests
1000X DTT Solution	1 x 100µl
1000X Protease Inhibitor Cocktail (PIC)	1 x 100µl
10X Pre-Extraction Buffer	1 x 10ml
Extraction Buffer	1 x 10ml
Extraction Cleaner	1 x 100µl
Extraction Pre-Cleaner	1 x 1ml

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.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