

ENE2 (Extraction Buffer) ab203377

Overview

Product name	ENE2 (Extraction Buffer)
Product overview	ENE2 (Extraction Buffer) (ab203377) is a part of the Nuclear Extraction Kit (ab113474). Refer to the kit protocol booklet for instructions of use. Please contact our Scientific Support team for more lot specific information.

Properties

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

Components	10 ml
ENE2 (Extraction Buffer)	1 x 10ml

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