

## Product datasheet

# Whole Cell Extraction Kit ab113475

[4 References](#) [1 Image](#)

### Overview

<b>Product name</b>	Whole Cell Extraction Kit
<b>Sample type</b>	Adherent cells, Suspension cells
<b>Assay time</b>	0h 45m
<b>Species reactivity</b>	<b>Reacts with:</b> Mouse, Human

**Predicted to work with:** Mammals 

### Product overview

Abcam's Whole Cell Extraction Kit (ab113475) provides a simple and selective method for extracting whole cell proteins in just 45 minutes for a variety of applications such as western blotting, cellular enzyme assays, or any other procedures requiring cellular proteins. ab113475 can be used to extract cellular proteins from mammalian cells and tissue samples.

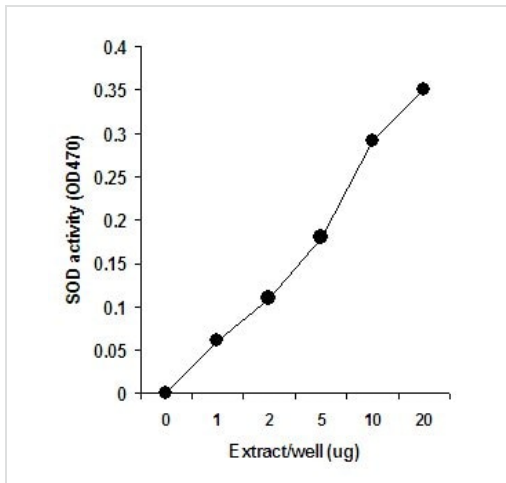
Not sure if this is the right product for you? Check out our [Sample Preparation Guide](#) for help.

### Properties

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

Components	100 tests
1000X DTT Solution	1 x 100µl
1000X Protease Inhibitor Cocktail	1 x 100µl
5X Extraction Buffer	1 x 20ml

### Images



Whole cell extracts were prepared from MCF-7 cells using ab113475. Activity of SOD was measured using different amounts of the extract.

Functional Studies - Whole Cell Extraction Kit  
(ab113475)

**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