

Product datasheet

Mammalian Cell Lysis Buffer 5X ab179835

17 References

Overview

Product name	Mammalian Cell Lysis Buffer 5X
Sample type	Tissue, Adherent cells, Suspension cells
Product overview	Mammalian Cell Lysis Buffer 5X (ab179835) is widely used to prepare mammalian cell and tissue lysates for use in a variety of downstream biochemical assays, especially those for quantification of enzymatic activity. This buffer does not contain SDS, a reagent that interferes with many activity assays.

The lysis buffer has been optimized for mammalian samples, but it can also be used for preparation of plant or bacterial samples.

The buffer just requires a simple 5-fold dilution, and minimal hands-on-time. Cell lysates are ready in only 15 minutes.

Notes	Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances. It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.
--------------	---

Properties

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

Components	10 ml
Mammalian Cell lysis buffer 5X	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