

Mammalian Cell Lysis Buffer 5X ab179835

17 References

製品の概要

製品名	Mammalian Cell Lysis Buffer 5X
サンプルの種類	Tissue, Adherent cells, Suspension cells
製品の概要	<p>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.</p> <p>The lysis buffer has been optimized for mammalian samples, but it can also be used for preparation of plant or bacterial samples.</p> <p>The buffer just requires a simple 5-fold dilution, and minimal hands-on-time. Cell lysates are ready in only 15 minutes.</p>
特記事項	<p>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.</p> <p>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.</p>

製品の特性

保存方法	Store at +4°C. Please refer to protocols.
------	---

内容	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.co.jp/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors