



Product datasheet

Mammalian Cell Extraction Kit ab65399

★★★★★ 1 Abreviews 19 References

概述

|      |  |
|------|--|
| 产品名称 | Mammalian Cell Extraction试剂盒   |
| 样品类型 | Tissue, Adherent cells, Suspension cells   |
| 检测类型 | Direct   |
| 检测时间 | 0h 20m   |
| 产品概述 | Abcam's Mammalian Cell Extraction Kit provides optimized cell extraction buffer, protease inhibitor cocktail, and DTT for convenient extraction of mammalian proteins from cultured cells and tissue samples, under nondenaturing conditions. Cell lysate prepared using the kit can be used in a variety of applications, such as enzyme activity assays (e.g. caspase activity assays), Western blot analysis, and others. The entire procedure takes less than 20 minutes.<br>Visit our <a href="#">FAQs page</a> for tips and troubleshooting. |
| 说明   | This product is manufactured by BioVision, an Abcam company and was previously called K269 Mammalian Cell Extraction Kit. K269-500 is the same size as the 500 test size of ab65399.   |

性能

存放说明 Please refer to protocols.

| 组件                          | 500 tests |
|-----------------------------|-----------|
| Cell Extraction Buffer      | 1 x 100ml |
| DTT                         | 1 x 110µl |
| Protease Inhibitor Cocktail | 1 vial    |

相关性 Mammalian Cell Extraction methods allow extraction of mammalian proteins from cultured cells and tissue samples.

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