

Genomic DNA Isolation Kit ab65358

10 References

概述

产品名称	基因组DNA分离试剂盒
样品类型	Tissue, Adherent cells, Suspension cells
检测类型	Direct
检测时间	2h 00m
产品概述	<p>Abcam's Genomic DNA Isolation Kit provides a simple and convenient procedure for rapid isolation of genomic DNA from mammalian cells and tissue samples with high yield and purity. The novel method requires less than 90 minutes to prepare highly pure genomic DNA. The extracted genomic DNA is free from protein and RNA, and suitable for a variety of applications such as PCR, DNA hybridization, enzyme manipulation, cloning, Southern blot, and array-based experiments.</p> <p>Visit our FAQs page for tips and troubleshooting.</p>
说明	<p>This product is manufactured by BioVision, an Abcam company and was previously called K281 Genomic DNA Isolation Kit. K281-50 is the same size as the 50 test size of ab65358.</p>

性能

存放说明	Store at -20°C. Please refer to protocols.
------	--

组件	50 tests
Cell Lysis Buffer	2 x 1.8ml
Enzyme Mix (Lyophilized)	1 vial
TE Buffer I	1 x 1.5ml

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