

Genomic DNA Isolation Kit ab65358

10 References

製品の概要

製品名	Genomic DNA Isolation Kit
サンプルの種類	Tissue, Adherent cells, Suspension cells
アッセイタイプ	Direct
全工程の試験時間	2h 00m
製品の概要	<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>
特記事項	<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>

製品の特性

保存方法 Store at -20°C. Please refer to protocols.

内容	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.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