

Product datasheet

Genomic DNA Isolation Kit ab65358

3 References

Overview

Product name	Genomic DNA Isolation Kit
Sample type	Tissue, Adherent cells, Suspension cells
Assay type	Direct
Assay time	2h 00m
Product overview	<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>
Tested applications	Suitable for: Purification

Properties

Storage instructions Store at -20°C. Please refer to protocols.

Components	50 tests
Cell Lysis Buffer	2 x 2ml
TdT Enzymes	1 vial
TE Buffer	1 x 1.5ml

Relevance Genomic DNA is the full complement of nuclear DNA contained in the cell or organism genome.

Applications

Our [Abpromise guarantee](#) covers the use of **ab65358** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
Purification		Use at an assay dependent concentration.

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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