

Product datasheet

Genomic DNA Extraction Kit ab156900

[16 References](#) [1 Image](#)

Overview

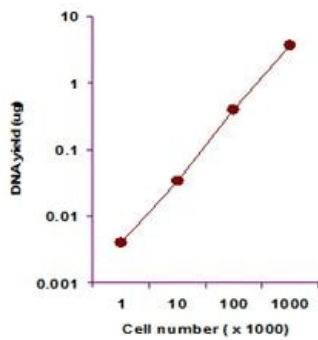
Product name	Genomic DNA Extraction Kit
Sample type	Adherent cells, Suspension cells, Whole Blood
Range	1 ng - 4000 ng
Assay time	0h 20m
Species reactivity	Reacts with: Mammals
Product overview	Genomic DNA Extraction Kit (ab156900) is designed for rapid isolation of pure genomic DNA from blood leukocytes or cultured mammalian cells in a small amount. The extracted DNA can be used for any molecular biology procedures such as PCR, restriction digestion, cloning and sequencing, etc. DNA yield can be up to 4 µg from 10 ⁶ blood leukocytes or cultured mammalian cells.

Properties

Storage instructions Please refer to protocols.

Components	50 tests	100 tests
DNG1 (Suspending Buffer)	1 x 16ml	2 x 16ml
DNG2 (DNA Digestion Solution)	1 x 1.1ml	1 x 2.2ml
DNG3 (DNA Digestion Powder)	1 vial	2 vials
DNG4 (DNA Capture Buffer)	1 x 16ml	2 x 16ml
DNG5 (DNA Elution Solution)	1 x 1ml	1 x 2ml
F-Collection Tube	50 units	100 units
F-Spin Column	50 units	100 units

Images



Genomic DNA was isolated from MCF7 cell line using ab156900.
The isolated DNA yield was quantified by real time PCR.

Genomic DNA purification using Genomic DNA
Extraction Kit (ab156900)

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.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