

Product datasheet

DNA Isolation Kit - Urine ab156899

[1 Image](#)

Overview

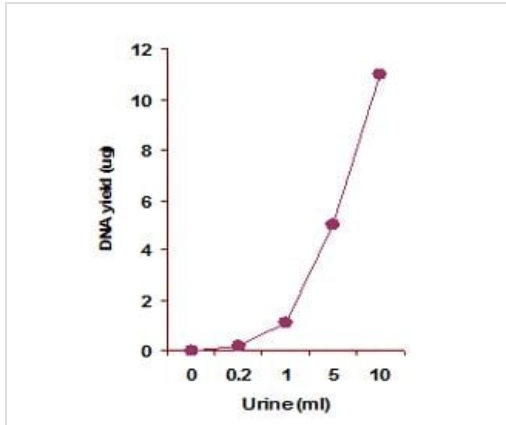
Product name	DNA Isolation Kit - Urine
Sample type	Urine
Range	1 ng - 2000 ng
Assay time	0h 20m
Product overview	The DNA Isolation Kit-Urine (ab156899) is a complete set of essential components which enables the experimenter to efficiently isolate DNA from urine. The entire procedure can be completed within 20 minutes and produces far superior results than any competitor kits.

Properties

Storage instructions Please refer to protocols.

Components	50 tests	100 tests
DNU1 (Suspending Buffer)	1 x 16ml	2 x 16ml
DNU2 (DNA Digestion Solution)	1 x 1.1ml	1 x 2.2ml
DNU3 (DNA Digestion Powder)	1 vial	2 vials
DNU4 (DNA Capture Buffer)	1 x 16ml	2 x 16ml
DNU5 (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 isolated from Human urine

Genomic DNA was isolated from Human urine using ab156899.
The isolated DNA yield was quantified by real time PCR.

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