

DNA Concentrator Kit ab156895

1 References

Overview

Product name	DNA Concentrator Kit
Sample type	DNA
Sensitivity	> 1 ng
Range	1 ng - 2000 ng
Assay time	0h 2m
Product overview	<p>The DNA Concentrator Kit (ab156895) is a complete set of essential components which enables the experimenter to efficiently concentrate DNA from various DNA samples with low concentration of DNA including those from microdissection samples and paraffin-embedded tissues. The entire procedure can be completed within 2 minutes and produces far superior results than any competitor kits.</p>

The kit is suitable for concentrating DNA from various DNA samples, especially from those with low concentration of DNA such as microdissection samples, paraffin-embedded tissues, restriction enzyme digestions, PCR reactions, and home-made minipreps.

Notes	<p>Concentrated DNA is highly pure and ready for direct use in PCR or other down- stream DNA manipulations such as ligation/transformation, sequencing, restriction digestion, microarray analysis, labeling, microinjection, and in vitro transcription.</p>
--------------	---

Properties

Storage instructions	Store at room temperature. Please refer to protocols.
-----------------------------	---

Components	50 tests	100 tests
DNC1 (Binding Buffer)	1 x 30ml	2 x 30ml
DNC2 (Wash Buffer)	1 x 2.5ml	2 x 2.5ml
DNC3 (DNA Elution Buffer)	1 x 1ml	1 x 2ml
F-Collection Tube	50 units	100 units

Components	50 tests	100 tests
F-Spin Column	50 units	100 units

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