

**ab156895**

# **DNA Concentrator Kit**

## **Instructions for Use**

For concentrating DNA from various low concentration DNA samples.

This product is for research use only and is not intended for diagnostic use.

# Table of Contents

---

Table of Contents	1
1. Background	2
2. Assay Summary	3
3. Materials Supplied	3
4. Storage and Stability	4
5. Materials Required, Not Supplied	4
6. Reagent Preparation	4
7. Assay Procedure	4

# 1. Background

---

The DNA Concentrator Kit uses a unique procedure and composition to efficiently concentrate DNA from various DNA samples with low concentration of DNA including those from microdissection samples and paraffin-embedded tissues. The kit has the following features:

- The fastest procedure available (< 2 minutes).
- Concentrated DNA is ultra-pure with efficient removal of contaminants (salts, proteins, enzymes, nucleotides, and primers).
- User-friendly binding solution and without chaotropic or other irritant reagents.

Abcam's DNA Concentrator 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.

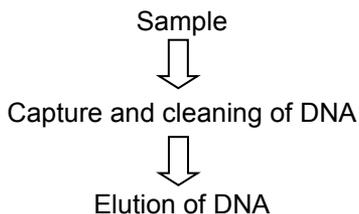
DNA size is from 70 bp to 20 kb; DNA quantity from 1 ng to 2 µg, optimal between 10 ng and 1 µg.

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.

ab156895 simply applies our proprietary DNA binding buffer to DNA sample. After transferring the sample to the specially designed Fast-Spin Column and washing it with the washing buffer, the DNA is easily recovered in 8-10 µL with elution solution. Concentrated DNA is ready for down-stream application.

## 2. Assay Summary

---



## 3. Materials Supplied

---

Item	50 tests	100 tests	Storage (Before Preparation)
DNC1 (Binding Buffer)	30 mL	2 x 30 mL	RT
DNC2 (Wash Buffer)	2.5 mL	2 x 2.5 mL	RT
DNC3 (DNA Elution Buffer)	1 mL	2 mL	RT
F-Spin Column	50	100	RT
F-Collection Tube	50	100	RT

## 4. Storage and Stability

---

The DNA Concentrator kit can be stored at room temperature (15-22°C).

## 5. Materials Required, Not Supplied

---

- Desktop centrifuge (up to 12,000 rpm)
- Pipettes and pipette tips
- 1.5 mL microcentrifuge tubes
- Ethanol (100%)

## 6. Reagent Preparation

---

Prepare final wash buffer:

Before starting, prepare the final wash buffer by adding 20 mL of 100% ethanol to DNC2.

## 7. Assay Procedure

---

**Note:** Always close spin columns before placing them in the microcentrifuge.

- a) Add 2 volumes of DNC1 to each volume of the DNA sample, and mix well. For example, add 200  $\mu$ L of DNC1 to 100  $\mu$ L of DNA sample. If required volume is less than 100  $\mu$ L, add 100  $\mu$ L of DNC1.
- b) Transfer the mixture to the column. Centrifuge at 12,000 rpm for 15 seconds. Discard the flowthrough. Replace the column to the collection tube

**Note:** *Maximum volume of the column is 600  $\mu$ L.*

- c) Add 200  $\mu\text{L}$  of DNC2 (final wash buffer) to the column and centrifuge at 12,000 rpm for 15 seconds. Discard the flowthrough and replace the column to the collection tube.
- d) Add 200  $\mu\text{L}$  of DNC2 to the column and centrifuge at 12,000 rpm for 30 seconds.
- e) Place the column in a new 1.5 mL vial. Add 6-8  $\mu\text{L}$  of DNC3 directly to the column filter, centrifuge at 12,000 rpm for 20 seconds to elute DNA. DNA is now ready for use or storage at  $-20^{\circ}\text{C}$ .



**UK, EU and ROW**

Email: [technical@abcam.com](mailto:technical@abcam.com) | Tel: +44-(0)1223-696000

**Austria**

Email: [wissenschaftlicherdienst@abcam.com](mailto:wissenschaftlicherdienst@abcam.com) | Tel: 019-288-259

**France**

Email: [supportscientifique@abcam.com](mailto:supportscientifique@abcam.com) | Tel: 01-46-94-62-96

**Germany**

Email: [wissenschaftlicherdienst@abcam.com](mailto:wissenschaftlicherdienst@abcam.com) | Tel: 030-896-779-154

**Spain**

Email: [soportecientifico@abcam.com](mailto:soportecientifico@abcam.com) | Tel: 911-146-554

**Switzerland**

Email: [technical@abcam.com](mailto:technical@abcam.com)

Tel (Deutsch): 0435-016-424 | Tel (Français): 0615-000-530

**US and Latin America**

Email: [us.technical@abcam.com](mailto:us.technical@abcam.com) | Tel: 888-77-ABCAM (22226)

**Canada**

Email: [ca.technical@abcam.com](mailto:ca.technical@abcam.com) | Tel: 877-749-8807

**China and Asia Pacific**

Email: [hk.technical@abcam.com](mailto:hk.technical@abcam.com) | Tel: 108008523689 (中國聯通)

**Japan**

Email: [technical@abcam.co.jp](mailto:technical@abcam.co.jp) | Tel: +81-(0)3-6231-0940

[www.abcam.com](http://www.abcam.com) | [www.abcam.cn](http://www.abcam.cn) | [www.abcam.co.jp](http://www.abcam.co.jp)