

AMPIXTRACT® Blood and Cultured Cell DNA Extraction Kit

Suitable for isolating DNA from blood leukocytes and cultured mammalian cells

The AMPIXTRACT® Blood and Cultured Cell DNA Extraction Kit is a complete set of essential components designed for rapid isolation of pure genomic DNA. The researcher can quickly and efficiently isolate DNA from blood leukocytes and cultured mammalian cells. The entire procedure can be completed within 20 minutes.

Ordering Information

[Order Online »](#)

ENZ-GEN503-0050	50Reactions
-----------------	-------------

Manuals, SDS & CofA

[View Online »](#)

- **Fast** - Entire procedure can be finished in 20 minutes with consistent isolation conditions.
- **Effective** - High efficiency of DNA isolation from blood leukocytes or cultured mammalian cells.
- **Reliable** - Our specially designed F-Spin Columns allow for DNA to be conveniently recovered.
- **Safe** - Use of non-toxic reagents and no phenol chloroform.

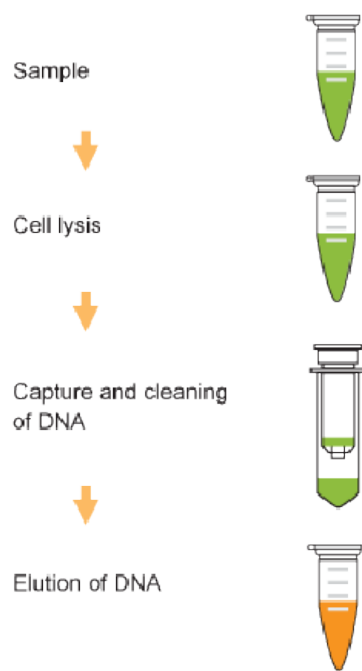


Figure 1. Schematic Procedure for using the AMPIXTRACT® Blood and Cultured Cell DNA Extraction Kit

Handling & Storage

Use/Stability The kit can be stored at room temperature (15-22°C) for up to 6 months from the date of shipment, with the exception of CB3. Upon receipt, CB3 should be stored at -20°C, or stored at 4°C as soon as it is dissolved in CB2 (stable for up to 6 months).

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Application Notes For isolating DNA from blood leukocytes and cultured mammalian cells

Assay Time 20 minutes

Contents

- CB1 (Suspending Buffer)
- CB2 (DNA Digestion Solution)
- CB3 (DNA Digestion Powder)
- CB4 (DNA Capture Buffer)
- CB5 (DNA Elution Solution)
- F-Spin Column
- F-Collection Tube
- Product Manual

Technical Info / Product Notes Input amount of blood leukocytes or cultured mammalian cells should be between 1 ng to 4 µg, with the optimal range between 10 ng and 1 µg. This kit can easily yield 1 ng to 4 µg of DNA in spin column format, which can be used for any molecular biology procedures such as PCR, restriction digestion, cloning and sequencing, etc.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)