

AMPIXTRACT[®] DNA Isolation Kit for Plasma/Serum

**Fastest procedure available to safely isolate
DNA from plasma/serum.**

AMPIXTRACT[®] DNA Isolation Kit for Plasma/Serum provides a simple method to isolate DNA from plasma, serum, and body fluids.

Ordering Information

[Order Online »](#)

ENZ-45018-0050	50Reactions
----------------	-------------

Manuals, SDS & CofA

[View Online »](#)

- Rapid - results in less than 15 minutes
- Safe - uses non-toxic reagents and no phenol chloroform
- High Efficiency - isolate DNA from plasma or serum containing as low as 0.1ng DNA

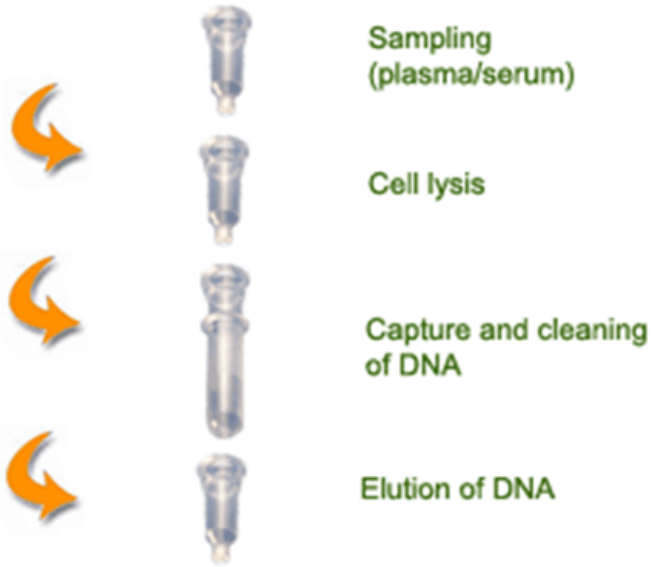


Figure 1. Schematic Procedure for using the AMPIXTRACT[®] DNA Isolation Kit for Plasma/Serum

Handling & Storage

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Application Notes For the isolation of DNA from plasma, serum, and body fluids.

Assay Time <15 min

Contents Digestion Solution, Digestion Powder, DNA Isolation Buffer, DNA Elution Solution, Spin Column, Collection Tube

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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