

AMPIXTRACT™

Magnetic Bead DNA Extraction Kit

Magnetic bead-based DNA extraction kit optimized for residual host cell DNA recovery

The AMPIXTRACT™ Magnetic Bead DNA Extraction Kit provides an efficient magnetic bead-based method for extracting residual host cell DNA from biopharmaceutical samples. Designed as the front-end sample preparation step for the AMPIGENE® Host Cell DNA Detection Kits (qPCR and dPCR), this kit delivers high-purity DNA suitable for sensitive downstream quantification. The streamlined protocol uses Proteinase K digestion followed by magnetic bead capture to efficiently recover trace-level DNA from complex matrices. The kit is compatible with all our host cell DNA detection workflows.

Residual host cell DNA (HCD) is a critical process-related impurity in biopharmaceutical manufacturing. Regulatory agencies including the FDA, EMA, and WHO require quantitative monitoring of residual HCD to ensure patient safety. Efficient DNA extraction from complex biopharmaceutical matrices is essential for accurate downstream detection by qPCR or dPCR methods. The magnetic bead-based approach provides consistent recovery across sample types while minimizing PCR inhibitor carryover.

Ordering Information

[Order Online »](#)

ENZ-GEN514-0100	100 extractions - 100extractions
-----------------	----------------------------------

Manuals, SDS & CofA

[View Online »](#)

- Magnetic bead-based extraction for consistent DNA recovery
- Optimized for biopharmaceutical matrices (mAbs, recombinant proteins)
- Compatible with both qPCR and dPCR downstream detection
- Works with CHO and E. coli host cell DNA workflows
- Proteinase K digestion for efficient protein removal
- Streamlined protocol — results in under 2 hours

Handling & Storage

Use/Stability	When stored and handled as directed, the kit is stable until the expiration date on the label.
Handling	Store Binding Buffer, Wash Buffer, and Elution Buffer at room temperature and rest of the components stored at 4°C. Add isopropanol to Binding Buffer and ethanol to Wash Buffer before first use per manual instructions.
Short Term Storage	4 °C
Long Term Storage	See component list
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application	DNA Extraction
Contents	Binding Buffer (40 mL); Wash Buffer (40 mL); 5X Proteinase K Buffer (3 mL); Proteinase K 20 mg/mL (1 mL); Magnetic Beads (5 mL); Elution Buffer (5 mL)
Technical Info / Product Notes	For use with AMPIGENE® CHO DNA Detection Kit (qPCR: ENZ-GEN516, dPCR: ENZ-GEN515) and AMPIGENE® E. coli DNA Detection Kit (qPCR: ENZ-GEN518, dPCR: ENZ-GEN517). See individual detection kit manuals for platform-specific extraction protocol modifications.



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)