

Automated nucleic acid extraction methods - Magnetic bead-based

	Method	Kit vendor	Fluent®	DreamPrep® NAP	Freedom EVO®	Resolvex® A200	Sample type
DNA Extraction	ZymoBIOMICS™ 96 MagBead DNA	Zymo Research®		○			Feces, soil, water, biofilms, swabs, saliva, body fluids (microbiomics/metagenomics)
	Quick-DNA™ MagBead Plus	Zymo Research		○			Biological fluids, blood, saliva, solid tissues, swabs and more
	NucleoMag® 96 Plant	Macherey-Nagel®			○		Plants
	Maxwell® HT 96 gDNA Blood Isolation System	Promega®	○				Blood, saliva
	ReliaPrep® Blood gDNA Miniprep System	Promega			○		Cells and tissue samples
	Wizard® SV 96 Genomic DNA Purification System	Promega			○		Rodent tail, cultured cells, human/animal cells and tissues
	Mag-Bind® Blood & Tissue DNA HDQ 96	Omega Bio-tek®	○				Whole blood, tissue, saliva, buccal swabs, cultured cells, mouse tail snips
	Mag-Bind Universal Pathogen	Omega Bio-tek	○				Tissue, urine, serum, and fecal samples
Cell-free DNA	Mag-Bind cfDNA	Omega Bio-tek	○			Serum, plasma	
Fragmented DNA	Mag-Bind TotalPure NGS	Omega Bio-tek	○			PCR products & NGS	
Plasmid DNA	Wizard MagneSil Tfx™ System	Promega			△		Bacterial culture
RNA isolation	Direct-zol™-96 MagBead RNA	Zymo Research				○	Cells, tissues, serum, plasma, blood, biological liquids, etc.

Legend: ○ Method qualified/verified by vendor¹ △ Method qualified/verified by Tecan¹ ◇ Method custom developed²



Method	Kit vendor	Fluent®	DreamPrep® NAP	Freedom EVO®	Resolvex® A200	Sample type
Viral DNA/RNA	NucleoMag Pathogen	Macherey-Nagel			○	Clinical samples
	Mag-Bind Viral DNA/RNA 96	Omega Bio-tek	○	○		Whole blood, serum, plasma, bodily fluids
	Maxwell® HT Viral TNA, Custom	Promega	◇		◇	Universal transport medium, serum, plasma and others
	QIAprep&® Viral RNA UM Kit	Qiagen	○			Swabs, saliva
	MagMax™ Viral/Pathogen Nucleic Acid Isolation	Thermo Fisher Scientific	◇		◇	Blood, swabs, urine, and viral transport media
	Quick-DNA/RNA™ Viral MagBead	Zymo Research	○	○	◇	Plasma, serum, urine, cell culture media, blood, saliva, cellular suspension, swab, fecal sample

Legend: ○ Method qualified/verified by vendor¹ Δ Method qualified/verified by Tecan¹ ◇ Method custom developed²

¹ Automated method developed according to the kit vendor's published protocol. Nucleic acids prepared with these methods meet the kit vendor's specifications and/or perform comparably to those prepared manually. Performance of the extraction procedure and the quality of the extracted nucleic acids was assessed either by the vendor (vendor verified) and/or by Tecan (Tecan verified).

² Automated method developed according to specific customer need. Extraction procedure tested by the customer.



Automated nucleic acid extraction methods - Vacuum-based

	Method	Kit vendor	Fluent®	DreamPrep® NAP	Freedom EVO®	Resolvex® A200	Sample type
DNA Extraction	NucleoSpin® 96 Blood	Macherey-Nagel®			○		Blood, biological samples
	NucleoSpin Food	Macherey-Nagel			○		Food, feed
	NucleoSpin Plant	Macherey-Nagel			○		Plants
	NucleoSpin Tissue	Macherey-Nagel			○	○	Animal tissue, cells, human tissue
	NucleoSpin DNA RapidLyse	Macherey-Nagel			○	○	Various tissue samples (mouse hard liver, kidney)
Plasmid DNA	NucleoSpin Plasmid	Macherey-Nagel			○		Bacteria, <i>E. coli</i>
	ChargeSwitch® Plasmid	Thermo Fischer Scientific®			△		<i>E. coli</i>
	ChargeSwitch-Pro Plasmid Miniprep	Thermo Fischer Scientific			△		Bacterial culture
	QIAprep® 96 Plus	Qiagen®				○	<i>E. coli</i>
	QIAprep 96 Turbo	Qiagen				○	<i>E. coli</i> , <i>S. cerevisiae</i> , <i>B. subtilis</i> , <i>A. tumefaciens</i>
RNA isolation	QIAamp® Viral RNA	Qiagen				○	Cell-free body fluids
	PureLink™ Pro 96 total RNA Purification	Thermo Fisher Scientific			○		Bacteria, yeast, plant, mammalian samples

Legend: ○ Method qualified/verified by vendor¹ △ Method qualified/verified by Tecan¹ ◇ Method custom developed²



Method	Kit vendor	Fluent®	DreamPrep® NAP	Freedom EVO®	Resolvex® A200	Sample type
Viral DNA/RNA	NucleoSpin RNA Virus	Macherey-Nagel			○	Cell-free biological fluid, plasma, serum
	NucleoSpin RNA Virus	Macherey-Nagel			○	Cell-free biological fluid, plasma, serum
	QIAprep&® Viral RNA UM Kit	Qiagen	○			Swabs, saliva
	QIAamp Viral RNA	Qiagen	○		○	Cell-free body fluids
	Quick-RNA™ Viral 96	Zymo Research				○

Legend: ○ Method qualified/verified by vendor¹ △ Method qualified/verified by Tecan¹ ◇ Method custom developed²

¹ Automated method developed according to the kit vendor's published protocol. Nucleic acids prepared with these methods meet the kit vendor's specifications and/or perform comparably to those prepared manually. Performance of the extraction procedure and the quality of the extracted nucleic acids was assessed either by the vendor (vendor verified) and/or by Tecan (Tecan verified).

² Automated method developed according to specific customer need. Extraction procedure tested by the customer.



www.tecan.com

DISCLAIMER:

For Research Use Only – not for use in diagnostic procedures. Example methods have not been validated. Please make a copy of this method in order to customize it to accommodate and validate your workflow according to your desired intended use and laboratory protocols. Tecan makes no claims regarding the performance of the example methods.

Tecan Group Ltd. makes every effort to include accurate and up-to-date information within this publication; however, it is possible that omissions or errors might have occurred. Tecan Group Ltd. cannot, therefore, make any representations or warranties, expressed or implied, as to the accuracy or completeness of the information provided in this publication. Changes in this publication can be made at any time without notice. For technical details and detailed procedures of the specifications provided in this document please contact your Tecan representative. This publication may contain reference to applications and products which are not available in all markets. Please check with your local sales representative.

Tecan, Freedom EVO, Fluent, and DreamPrep are registered trademarks of Tecan Trading Ltd., Männedorf, Switzerland; Resolvex is a registered trademark of Tecan SP, Inc.; Maxwell, Magnesil, Promega and Wizard are registered trademarks of Promega Corp.; ChargeSwitch, PureLink and MagMax are registered trademarks of Life Technologies Corp.; Omega Bio-Tek is a registered trademark of Qi Guo, Mag-Bind is a registered trademark of Omega Bio-tek Inc.; NucleoMag and NucleoSpin are registered trademarks of Macherey, Nagel GmbH & Co.; QIAprep, QIAprep& and QIAamp are registered trademarks of Qiagen GmbH; Zymo Research, ZymoBIOMICS are registered trademarks of Zymo Research Corporation.

All mentioned trademarks are protected by law. In general, the trademarks and designs referenced herein are trademarks, or registered trademarks, of Tecan Trading Ltd., Männedorf, Switzerland. A complete list may be found at www.tecan.com/trademarks. Product names and company names that are not contained in the list but are noted herein may be the trademarks of their respective owners.

Australia +61 3 9647 4100 **Austria** +43 62 46 89 33 **Belgium** +32 15 42 13 19 **China** +86 21 220 63 206 **Denmark** +45 70 23 44 50 **France** +33 4 72 76 04 80 **Germany** +49 79 51 94 170 **Italy** +39 02 92 44 790 **Japan** +81 44 556 73 11 **Netherlands** +31 18 34 48 17 4 **Singapore** +65 644 41 886 **Spain** +34 93 490 01 74 **Sweden** +46 8 750 39 40 **Switzerland** +41 44 922 89 22 **UK** +44 118 9300 300 **USA** +1 919 361 5200 **Other countries** +43 62 46 89 33

