

Freedom EVO	Vendor qualified/verified	Tecan qualified
ChargeSwitch® Plasmid kit		●
Dynabeads® mRNA Purification kit		●
DNA IQ™ System and Casework Extraction	●	
Extract-N-Amp™ Plant Kit		●
IPureLink™ Pro 96 Total RNA Purification Kit		●
Maxwell® HT DNA Blood Kit	●	
Maxwell HT Viral tDNA Kit (custom)	●	
Maxwell HT Simply RNA kit (Paxgene, custom)	●	
NucleoSpin® 96 Blood Kit	●	
NucleoSpin 8/96 Food Kit	●	
NucleoSpin 96 Plant Kit	●	
NucleoSpin 8/96 Plasmid Kit	●	
NucleoSpin 96 RNA Kit	●	
NucleoSpin 96 Tissue Kit	●	
NucleoSpin 96 PCR Clean-up	●	
NucleoMag® Blood Kit (200 µl)	●	
NucleoMag Pathogen	●	
NucleoMag Plant Kit	●	
NucleoMag RNA Kit	●	
NucleoMag Tissue Kit	●	
NucleoMag Virus Kit	●	
NucleoMag VET	●	
Wizard® SV96 Genomic DNA Purification System	●	
Wizard® MagneSil® DNA Purification System/Tfx™	●	

Fluent	Vendor qualified/verified	Tecan qualified
Maxwell HT DNA Blood Kit	●	
Maxwell HT Viral tDNA Kit (custom)	●	
Maxwell HT Simply RNA kit (tissue, custom)	●	
NucleoSpin 96 Tissue Kit		●