



Choose the best solution for your environmental nucleic acid isolation

[Discover the microbiome workflow](#)



Isolating nucleic acids from environmental samples can be a challenging part of microbiome research. Our products help you isolate DNA and RNA from soil, water, air, biofilm, plants and seeds. The optimized bead beating technology in our kits ensures unbiased lysis of the toughest microbes. We include Inhibitor Removal Technology® to eliminate humic acids, polyphenols, polysaccharides, heavy metals and other impurities. This leaves you with the purest possible nucleic acids for downstream applications. To suit your processing needs, we offer kit throughputs from manual to 96-well for a variety of **automation platforms**.

Soil*

Kit	Starting amount	Recommendations for usage	Automated on	Cat. no.
DNA				
DNeasy® PowerSoil® Pro	250 mg	For all soil types	QIAcube® Connect (optional)	47014 47016
DNeasy PowerSoil Pro QIAcube	250 mg	For all soil types	QIAcube Connect	47126
DNeasy 96 PowerSoil Pro QIAcube HT**	250 mg	Up to 96 samples	QIAcube HT	47021
DNeasy 96 PowerSoil Pro	250 mg	Up to 96 samples	Manual processing on plates	47017
DNeasy PowerMax® Soil	Up to 10 g	Low biomass soil	Manual processing in 50 mL tubes	12988-10
MagAttract® PowerSoil Pro DNA**	100 mg	Up to 96 samples	KingFisher™ epMotion®	47109 coming soon
DNeasy PowerSoil Pro WGS SeqSet	250 mg	For all soil types	Automation compatible	47165, 47175, 47185, 47195, 47203
RNA				
RNeasy PowerMax Soil Pro	Up to 15 g		Manual processing in 50 mL tubes	47082

* For all soil types including rhizosphere, compost, clay, top soil, etc.

** Bead Plates/tubes must be purchased separately for QIAcube HT and MagAttract Kits: Cat No. 19301 PowerBead Pro Tubes (2 ml) (50) or Cat No. 19311 PowerBead Pro Plates (4). To learn how to properly use the bead plates, please watch our explainer video.

Water & Air

Kit	Starting sample	Recommendations for usage	Automated on	Cat. no.
DNA				
DNeasy PowerWater®	Water or Air Filter	For most filter membrane types	QIAcube Connect (optional)	14900-50-NF 14900-100-NF
DNeasy PowerWater Sterivex™	Sterivex Filter Unit	Sterivex Filter Units		14600-50-NF
RNA				
RNeasy PowerWater	Water or Air Filter	For most filter membrane types		14700-50-NF
DNA & RNA				
AllPrep® PowerViral® DNA/RNA	200 µl waste water	Viral nucleic acid isolation from waste water	QIAcube Connect (optional)	28000-50

Biofilm

Kit	Starting sample	Automated on	Cat. no.
DNA			
DNeasy PowerBiofilm®	50–200 mg	QIAcube Connect (optional)	24000-50
RNA			
RNeasy PowerBiofilm	50–200 mg		25000-50

Plants & Seeds

Kit	Starting sample	Recommendations for usage	Automated on	Cat. no.
DNA				
DNeasy Plant Pro	Up to 100 mg	For plant cells, tissues, seeds or fungi	QIAcube Connect (optional)	69204
DNeasy 96 Plant	10–50 mg	Up to 96 samples	Manual processing on plates	69181
RNA				
RNeasy PowerPlant	10–50 mg	For plants or seeds		13500-50



Visit www.qiagen.com/applications/microbiome or contact QIAGEN today to get more out of your environmental samples!

For up-to-date licensing information and product-specific disclaimers, see the respective QIAGEN kit handbook or user manual. QIAGEN kit handbooks and user manuals are available at www.qiagen.com or can be requested from QIAGEN Technical Services or your local distributor.

Trademarks: QIAGEN®, Sample to Insight®, QIAcube®, DNeasy®, RNeasy®, Inhibitor Removal Technology®, MagAttract®, PowerBiofilm®, PowerMax®, PowerPlant®, PowerSoil®, PowerViral®, PowerWater®, AllPrep®, (QIAGEN Group); KingFisher™ (Thermo Fisher Scientific, Inc.); EpMotion® (Eppendorf AG); Sterivex™ (Millipore). Registered names, trademarks, etc. used in this document, even when not specifically marked as such, may still be protected by law. QPRO-5156 09/ © 2023 QIAGEN, all rights reserved.

Ordering www.qiagen.com/shop | Technical Support www.support.qiagen.com | Website www.qiagen.com