



# Choose the best solution for your human microbiome samples

[Discover the microbiome workflow](#)



Isolating nucleic acids from human samples for microbiome research can be challenging. Our products help you isolate nucleic acids from various sample types, with flexible automation and throughput options. The integrated bead beating technology ensures unbiased lysis of the toughest microbes. The Inhibitor Removal Technology<sup>®</sup> eliminates PCR inhibitors such as lipids, heme, polyphenols and polysaccharides from fecal samples to ensure pure nucleic acids that work in downstream applications. To suit your processing needs, we offer kit throughputs from manual to 96-well for a variety of **automation platforms**.

Kit	Sample type					Automated on	Format	Cat. no.
	Fecal/gut	Biofluids	Swab	Biofilm	Waste water			
<b>Nucleic acid stabilization</b>								
PowerProtect DNA/RNA	●		●			—	Bulk solution	14810
<b>Microbial gDNA</b>								
QIAamp <sup>®</sup> PowerFecal <sup>®</sup> Pro DNA	●		●			QIAcube <sup>®</sup> Connect (optional)	Single column	51804
QIAamp PowerFecal Pro DNA QIAcube	●		●			QIAcube <sup>®</sup> Connect	Single column	51826
QIASymphony <sup>®</sup> PowerFecal Pro DNA	●		●			QIASymphony SP	Single column	938036
DNeasy <sup>®</sup> PowerBiofilm <sup>®</sup>				●		QIAcube Connect (optional)	Single column	24000-50
DNeasy 96 PowerSoil <sup>®</sup> Pro	●		●			Manual processing on plates	96-well plate	47017
DNeasy 96 PowerSoil Pro QIAcube HT*	●		●			QIAcube HT	96-well plate	47021
QIAamp DNA Microbiome**		●	●			QIAcube Connect (optional)	Single column	51704

\* 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.  
 \*\* Enables efficient depletion of human host gDNA

Kit	Sample type					Automated on	Format	Cat. no.
	Fecal/gut	Biofluids	Swab	Biofilm	Waste water			
<b>Microbial gDNA</b>								
MagAttract® PowerSoil Pro DNA*	●		●			KingFisher™ epMotion®	96-well plate	47109 47119
QIAamp PowerFecal Pro WGS SeqSet	●		●			Automation compatible	Single column + sequencing kit	51825, 51835, 51845, 51855, 51863
<b>Microbial RNA</b>								
RNeasy PowerFecal Pro	●		●			QIAcube Connect (optional)	Single column	78404
RNeasy PowerBiofilm				●			Single column	25000-50
<b>Microbial gDNA &amp; Microbial RNA</b>								
AllPrep PowerFecal Pro DNA/RNA	●		●				Single column	80254
AllPrep PowerViral® DNA/RNA	●				●	QIAcube Connect (optional)	Single column	28000-50
MagAttract PowerMicrobiome DNA/RNA	●		●			KingFisher™ epMotion®	96-well plate	27500-4-EP 27600-4-KF
<b>Host gDNA</b>								
QIAamp Fast DNA Stool Mini***	●		●			QIAcube Connect (optional)	Single column	51604

\* 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.  
 \*\*\* Captures host genomic DNA and easy-to-lyse microbial gDNA



Visit [www.qiagen.com/applications/microbiome](http://www.qiagen.com/applications/microbiome) or contact QIAGEN today to get more out of your fecal 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](http://www.qiagen.com) or can be requested from QIAGEN Technical Services or your local distributor.

Trademarks: QIAGEN®, Sample to Insight®, QIAamp®, QIAcube®, QIASymphony®, AllPrep®, DNeasy®, RNeasy®, Inhibitor Removal Technology®, MagAttract®, PowerFecal®, PowerBiofilm®, PowerMicrobiome®, PowerSoil®, PowerViral®, (QIAGEN Group); KingFisher™ (Thermo Fisher Scientific, Inc.) EpMotion® (Eppendorf AG). Registered names, trademarks, etc. used in this document, even when not specifically marked as such, may still be protected by law. QPRO-5169 2023© QIAGEN, all rights reserved.

Ordering [www.qiagen.com/shop](http://www.qiagen.com/shop) | Technical Support [www.support.qiagen.com](http://www.support.qiagen.com) | Website [www.qiagen.com](http://www.qiagen.com)