

Third-Party Reagent Protocols for Bravo NGS and NGS Workstation

The list below is a sample of readily available third-party protocols for Bravo. Custom Bravo protocols can be created for your specific needs. Contact your local field representative for more information.

Company	NGS Reagent	Bravo Options*
Illumina	COVIDSeq	A, B
	Illumina Standard Total RNA	A, B
	Illumina Standard mRNA	A, B
	Illumina DNA Prep with Enrichment	A, B
	Illumina DNA PCR-free Library Prep Tagmentation	A
	Nextera XT	A, B
	Nextera Flex	A, B
	Illumina Stranded Total RNA with RiboZero	A
Watchmaker	mRNA Library Preparation	B
KAPA Biosystems (Roche)	HTP Library Preparation Kit	A, B
	HyperPrep DNA Library Preparation Kit	A, B
	HyperPlus DNA Library Preparation Kit	A, B
	Stranded RNA-Seq Library Preparation Kit	B
	Stranded mRNA-Seq Library Preparation Kit	A, B
New England Biolabs (NEB)	NEBNext	A
	Low Input RNA	A
	Immune Sequencing Kit Human	A
	NEBNext DNA Library Prep	B
	NEBNext Ultra II RNA	A, B
Oxford Nanopore	General Ligation LSK114 workflow	B
Pacific Biosciences	SMRTbell Library Preparation (10 Kb) (PacBio RS II)	B
Clontech	Smart-Seqv4 Ultra Low RNA	A, B
10X Genomics	Chromium Next GEM Single Cell 5' v2	A
	Single Cell 3' V3,1 (Dual Index)	A
	Single Cell 3' V3,1	B

[*Bravo NGS \(Option A\) Bravo NGS Workstations \(Options B\)](#)