

Clinical Facilities

Vendor Qualified/Verified    Field Developed

Transfusion, Transplantation Medicine (HLA typing)	NGSgo		●
	TruSight® HLA v2.0 Sequencing Panel®	●	
Medical Microbiology, Virology, Hygiene, Infectious diseases	QiaSeq FX		●
	Nextera XT	●	
	NEBNext™ Ultra II DNA	●	
Oncology, Cancer Institute, Molecular Pathology	Oncomine Fucos Assay		●
	QIAseq™ Library Preparation		●
	TruSeq Custom Amplicon Low Input	●	
	Kapa™ HyperPlus		●
Human Genetics, Rare genetic diseases	SureSelect™ QXT Target Enrichment		●
	TruSeq RNA Access	●	
	NEBNext Ultra II RNA		●
	Nextera® Rapid Capture	●	
	NimbleGen™ SeqCap™ EZ Library Prep	●	
	Kapa HyperPrep		●
	Ion AmpliSeq™	●	