

Summary of Applications

A diverse array of applications can be automated using Automation Solutions platforms. The following table presents some examples of applications available. For other applications and for [professional services and consultation](#), contact [Automation Solutions](#).

Applications		Bravo Platform	BenchCel Workstation	BioCel System
Cell Biology	ADME-Tox	✓	✓	✓
	Cell Viability	✓	✓	✓
	Cell-Based Assays	✓	✓	✓
Genomics	Genotyping			✓
	DNA Isolation	✓	✓	✓
	Microarray	✓	✓	✓
	RNA Isolation	✓	✓	✓
	PCR Preparation/Purification	✓	✓	✓
Immunoassays	ELISA		✓	✓
Screening	Cell-Based Assays	✓	✓	✓
	Compound Management		✓	✓
	Enzyme Assays		✓	✓
	Microplate Replication and Reformatting	✓	✓	✓
	Sample Picking		✓	✓
Liquid Handling	Magnetic Bead Purification	✓	✓	✓
	Microplate Filtration	✓	✓	✓
	Serial Dilution	✓	✓	✓
	Tip Washing	✓	✓	✓
Microplate Management	Centrifugation		✓	✓
	Barcode Tracking	✓	✓	✓
	Labeling		✓	✓
	Transfer	✓	✓	✓
	Sealing		✓	✓
	Storage		✓	✓