

Compatible Microarray Scanners for Imaging scWest Chips

Scanner Name	Laser Count	Company
SpotLight Turbo	Up to 6	ArrayIT
SureScan	2 (simultaneous)	Agilent
InnoScan 710	2	Innopsys
InnoScan 910	2	Innopsys
InnoScan 1100AL	3	Innopsys
GenePix 4000B	2	Molecular Devices
GenePix 4100A	2	Molecular Devices
GenePix 4300A	Up to 4	Molecular Devices
GenePix 4400A	Up to 4	Molecular Devices
GenePix Professional 4200A	Up to 4	Molecular Devices
ScanArray Express	Up to 5	PerkinElmer
ScanArray Express HT	Up to 5	PerkinElmer
ScanArray 5000	4	PerkinElmer
LS Reloaded	Up to 4	Tecan
PowerScanner	2	Tecan
ArrayIT SpotLight	2	Telechem
G2565CA	2	Agilent

NOTE: Recommended scanners validated by ProteinSimple are shown in **bolded** text. Other scanners listed should be compatible based on specifications provided by manufacturer.

This list is not exhaustive so please ask your ProteinSimple representative if you have a scanner that's not listed here.