

Rev. Date November 9, 2015

Feature Extraction Compatibility Matrix for Non Agilent scanners

Manual gridding may be required for any 3rd party scanner.

Feature Extraction Standalone v12.0 and Feature Extraction for Cytogenomics v2.9 are able to support Agilent CGH array images generated on select non-Agilent scanners.

Scanner	CGH+SNP		CGH		GE		CHIP-CHIP/Methylation		miRNA	
	HD	G3	HD	G3	HD	G3	HD	G3	HD	G3
NimbleGen MS200/Tecan Power Scanner	Yes	Yes	Yes	Yes	Yes	Yes	Yes	Yes	Yes	Yes
Innopsys Innoscan 710 & 900	Yes	Yes	Yes	Yes						
Molecular Dynamics GenePix 4300 & 4400	Yes	Yes	Yes	Yes						

Note: No verification of algorithm performance or protocols was performed on non-CGH TIFF images for the Axon and Innopsys scanners.

The performance of the GenePix 4000 and 4100 scanners was not tested but they should perform similarly to the 4300/4400 models according to the manufacturer.

[Download Agilent CytoGenomics Software](#)

[Download Feature Extraction Standalone Software](#)

[Feature Extraction User Guide \(G4460-90053\)](#)

Recommended Scanner Settings

[Feature Extraction for CytoGenomics 2.5 - Settings for Roche NimbleGen MS200 Scanners \(G1662-90026\)](#)

[Feature Extraction for CytoGenomics 2.9 - Settings for Innopsys Innoscan 710 & 900 Scanners \(G1662-90036\)](#)

[Feature Extraction for CytoGenomics 2.9 - Settings for Molecular Dynamics GenePix 4300 & 4400 Scanners \(G1662-90035\)](#)

NimbleGen/TECAN, Innopsys, Molecular Devices Scanner Support

At this time, Agilent has performed feature extraction of aCGH scans from NimbleGen MS200, Innopsys Innoscan 710 and 900, and Molecular Devices GenePix 4300 and 4400 using Feature Extraction. Verification of algorithm performance or protocols have been performed.

© Agilent Technologies, Inc. 2014, 2015

G1662-90043 Revision C1, November 2015

For Research Use Only. Not for use in diagnostic procedures.