

Instrument	File type and # files exported
ABI PRISM [®] 7700	All data for all dyes are exported in 1 file (.tab)
ABI PRISM [®] 7900	All data for all dyes are exported in 1 file (.txt)
ABI PRISM [®] 7000, 7500	One file per curve type (Amplification, Melt), includes all dyes. (.csv)
Roche LightCycler [®] and LightCycler [®] 2.0	One file per curve type (amp, melt) per dye type (.txt)
BioRad iCycler [®] iQ	iCycler [®] files must be copied to Excel and saved as .txt files.* One file per curve type (amp, melt) per dye type (.txt)
Stratagene MX3000P [™]	All data for all dyes are exported in 1 file (.txt)
MJ Research DNA Engine Opticon [®] 2	All data for all dyes are exported in 1 file (.csv)
Cepheid SmartCycler [®] II	One file for all Amplification dyes, 1 file per dye for Melt data (.csv)

.tab or **.txt**: Tab-delimited file

.csv: Comma-separated File

* If the Plexor[™] Analysis Software is installed on the computer that is running the BioRad iCycler iQ[™] Software, data can be copied directly from the BioRad iCycler[™] software then pasted into the Plexor[™] File import screen. Select "BioRad iCycler iQ[™] (via paste)" as the instrument in the Assay Setup screen.