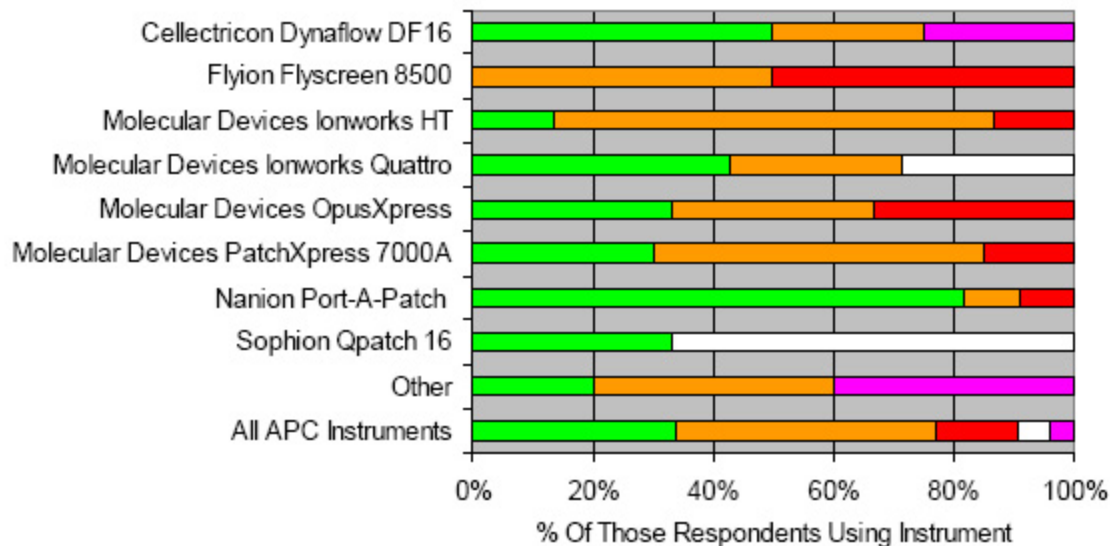


Respondent's Opinion on Overall APC Instrument Performance

Overall APC Instrument Performance



- High (very happy - exceeds expectations)
- Medium (acceptable)
- Low (not satisfied - needs improvement)
- Too Early To Say
- Can't Say

© HTStec 2005

Manufacturer/Model	Opinion On Overall APC Instrument Performance					No. Different Instrument Installs Assessed
	High (very happy - exceeds expectations)	Medium (acceptable)	Low (not satisfied - needs improvement)	Too Early To Say	Can't Say	
Cellectricon Dynaflo DF16	2	1	0	0	1	4
Flyion Flyscreen 8500	0	3	3	0	0	6
Molecular Devices Ionworks HT	2	11	2	0	0	15
Molecular Devices Ionworks Quattro	3	2	0	2	0	7
Molecular Devices OpusXpress	1	1	1	0	0	3
Molecular Devices PatchXpress 7000A	6	11	3	0	0	20
Nanon Port-A-Patch	9	1	1	0	0	11
Other	1	2	0	0	2	5
Sophion Qpatch 16	1	0	0	2	0	3
TOTAL	25	32	10	4	3	74