

SOFTWARE	INTERFACE	AMPLIFIERS
Clampex9	Digidata	all manual amplifiers* Multiclamp 700 A/B**
Pulse	ITC 16 ITC 18 LIH1600	all manual amplifiers* EPC 9/10
PatchMaster	ITC 16 ITC 18 LIH 1600	all manual amplifiers* EPC 9 EPC 10
GePulse	ITC 16 ITC 18 National Instruments	all manual amplifiers*

* e.g. Axopatch 200B, EPC 7, PC-ONE

** Software interface to MulticlampCommander is available