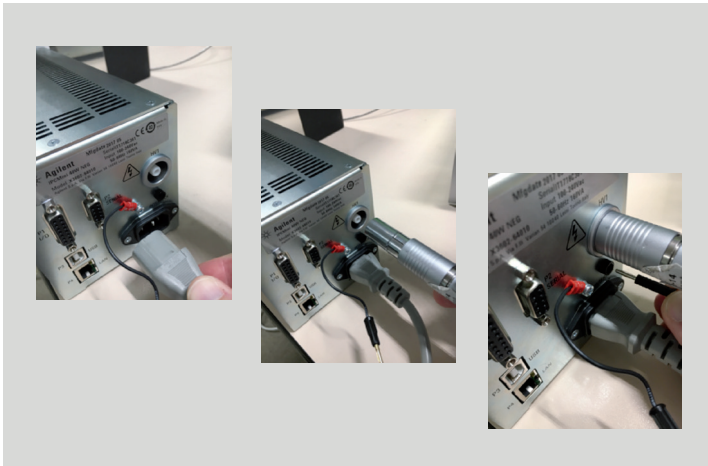


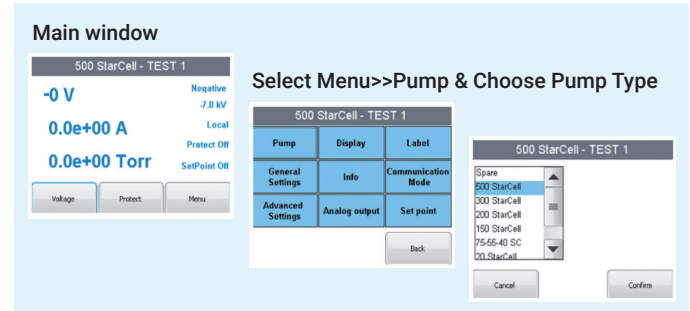
# Agilent IPCMini Quick Start Guide



## Connect HV and AC Power



## Select Pump Type



Press & Hold  
HV On/Off Switch



