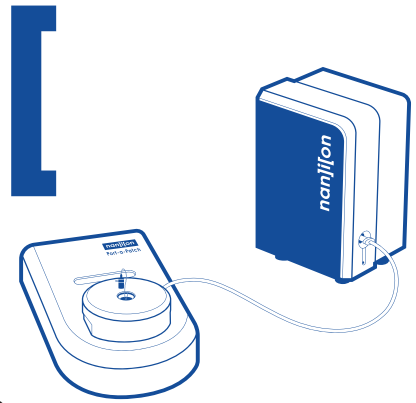


Port-a-Patch Configurations



Key Features

- 1 Easy-To-Use, Even For Non-Experts
- 2 Voltage & Current Clamp as Standard
- 3 Whole-Cell & Single Channel Recordings
- 4 External & Internal Perfusion
- 5 Temperature Control (Cooling & Heating)
- 6 Compatible With Most Amplifiers
- 7 Light Stimulation (Optional)
- 8 Full Control Over Experimental Parameters For Fast Assay Development

Packages	Premium	Perfusion	Basic	Light	Mini
Amplifier	HEKA EPC 10 Also Compatible With Axopatch 200B or MultiClamp 700A & B	HEKA EPC 10 Also Compatible With Axopatch 200B or MultiClamp 700A & B	HEKA EPC 10 Also Compatible With Axopatch 200B or MultiClamp 700A & B	No Amp. Included Also Compatible With HEKA EPC 10, Axopatch 200B, or MultiClamp 700A & B	Integrated Tecella Pico & Headstage
Features	Internal Perfusion External Perfusion Temperature Control Suction Control Pro Unit	External Perfusion Temp. Control Suction Control Unit	Suction Control Unit <i>Upgrades for Perfusion & Temp. Control Available</i>	Suction Control Unit <i>Upgrades for Perfusion & Temp. Control Available</i>	Suction Control Unit
Data Acquisition Software	Patchmaster pClamp Suite	Patchmaster pClamp Suite	Patchmaster pClamp Suite	Patchmaster pClamp Suite	WinWCP
Consumables	500 NPC-1 Chips Reagent Kit Maintenance Kit Electrode Service Kit Spare Part Kit	500 NPC-1 Chips Reagent Kit Maintenance Kit	250 NPC-1 Chips Reagent Kit Maintenance Kit	250 NPC-1 Chips Reagent Kit Maintenance Kit	50 NPC-1 Chips Reagent Kit
Warranty & Service Plan	1 Year Warranty 1 Year Service Plan	1 Year Warranty	1 Year Warranty	1 Year Warranty	1 Year Warranty
Computer & Monitor Included	✓	✓	✓	✓	Optional
Voltage Clamp	✓	✓	✓	✓	✓
Current Clamp	✓	✓	✓	Dependent on Integrated Amplifier	✓



Accelerate Your Research

The Port-a-Patch is a small and easy-to-use complete patch clamp setup with multiple add-ons available to make sure you get the right configuration for your application.