

Midi Plus Pipette Controller for Use in Biosafety Cabinet

The Company operates one of the world's largest industrial network of vaccine production with nearly 9,000 employees dedicated to vaccine development and exporting to 150+ countries.

Challenge

- The serological pipettors will be used in a barrier technology and must be designed as to allow easy decontamination
- Maximum 5h in discontinue use based on estimated time of sampling of 1 min (From 4 to 100 samples)
- Fixation can be added to support pipette controller location in the cabinet

The Solution

- Midi Plus can run ~2000 pipetting cycles when fully charged, depending on dispensing method, used liquid and volumes.
- Autoclavable nose cone, silicone adapters and 0.2 μm venting filter (121°C, 20 min, 1 bar/100 kPa). Can be cleaned from the outside with 70% ethanol or mild cleaning agents
- Integrated fold-out stand

The Value

- Optimized battery allows for continuous use during user work shift
- Easy cleaning to keep cross-contamination under control
- Fold-out stand saves bench space when Midi Plus is not in use and keeps the pipette free from the table's surface to avoid contamination



For more information,
visit www.sartorius.com

SARTORIUS