

Speedcal Mobile

Fast Calibration of Multi-Channel Pipettes



Product Information

The Speedcal Mobile system was developed to meet the specific requirements for calibrating multi-channel pipettes. Its innovative design and sophisticated technology simplify the entire calibration process. Parallel connection of up to 12 balances, which simultaneously determine the dispensed volumes from a multi-channel pipette, make Speedcal Mobile the fastest and most effective system on the market for this task. The integrated Web Service interface enables Speedcal Mobile to be connected to all pipette calibration software. Speedcal Mobile is suitable for long-term use in laboratories and, thanks to its practical carrying case, also for use in mobile service on site.

Benefits

- Testing twelve channels as per DIN EN ISO 8655 10.3 within approximately 10 minutes saves time
- Unique system, available with either 4, 8, or 12 channels, with the option of subsequently adding up to 12 channels
- Conforms to all national (ISO17025) and international (ISO8655) regulations



Technical Specifications

Overview of the Sartorius Multi-Channel Pipette Calibration System

Model	Speedcal-M12	Speedcal-M8	Speedcal-M4
Number of channels	12	8	4
Sartorius weigh cells	12	8	4
Weighing capacity	21 g per channel	21 g per channel	21 g per channel
Resolution	0.01 mg	0.01 mg	0.01 mg
Taring range	-21 g per channel	-21 g per channel	-21 g per channel
Repeatability Standard deviation	0.02 mg	0.02 mg	0.02 mg
Stabilization time	<4 s	<4 s	<4 s
Sensitivity to temperature drift	1 ppm/°K	1 ppm/°K	1 ppm/°K
Max. volumes	21 ml per channel	21 ml per channel	21 ml per channel
Interface	Web Service	Web Service	Web Service
Weighing module dimensions	450 × 270 × 150 mm	450 × 270 × 150 mm	450 × 270 × 150 mm
Weight of weighing module	14 kg	12 kg	10 kg
E-box dimensions	600 × 450 × 150 mm	600 × 450 × 150 mm	600 × 450 × 150 mm
E-box weight	5 kg	5 kg	5 kg
Case transport weight	~ 45 kg	~ 45 kg	~ 45 kg
Case transport dimensions	610 × 485 × 608 mm	610 × 485 × 608 mm	610 × 485 × 608 mm
Pallet transport weight	~ 95 kg	~ 93 kg	~ 90 kg
Pallet transport dimensions	800 × 665 × 910 mm	800 × 665 × 910 mm	800 × 665 × 910 mm




Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Straße 20
37079 Göttingen
Phone +49 551 308 0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906

 For further information, visit
www.sartorius.com