

Save Costs, Time and Calibrate More Pipettes with SpeedCal® Mobile



Calculate Your Savings, Compared to the Fastest Competitor Devices

Pipette Model	How many Multi-Channel Pipettes do you calibrate per day?	How fast is your current system?	Labor costs in \$/h	Working days per year
12-Channel				
8-Channel				
4-Channel				
SpeedCal® Mobile	Saved Labor Time per Year in Hours	Labor Costs Saved perYear	Up to x more 12-Channel Calibrations per Year	
M4				
M12				

All blue fields are editable fields. Once filled, they will show your saved labor time in hours per year, your saved laborcosts, and how many more 12 Channel Pipettes you could calibrate when switching to a SpeedCal®-M4 or M12.

What are Your Benefits?



Save on Labor Time

Free up annoying calibration time for other work in your lab



Increase Your Throughput and Grow Your Business

SpeedCal®'s high-speed multi-channel calibration lets you handle more pipettes per year, enabling higher turnover and expanding your operational capacity.



Work Smarter, at Lower cost

SpeedCal® simplifies the calibration workflow so your lab can operate more efficiently, reducing training time and supporting cost-effective daily operations.

Learn More:



SpeedCal® delivers ultra-fast, compliant multi-channel pipette calibration. It boosts throughput, reduces training effort, and adapts to your growth with upgradeable configurations. More pipettes calibrated. Less time wasted. Maximum efficiency for modern calibration labs.