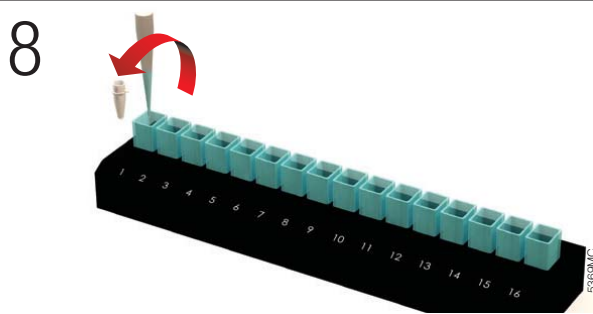
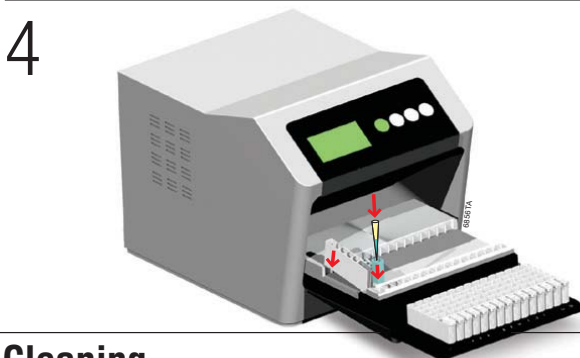
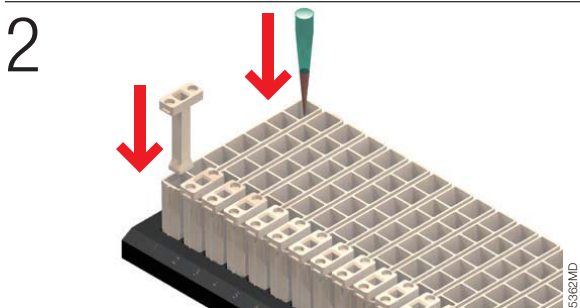
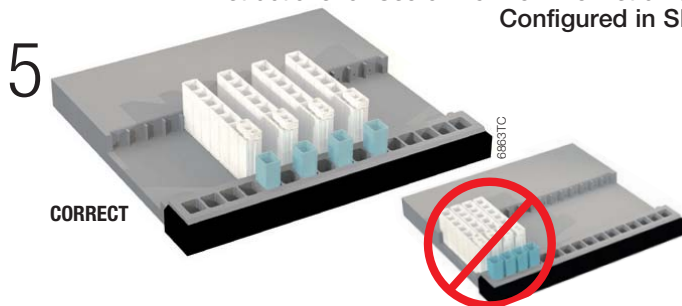
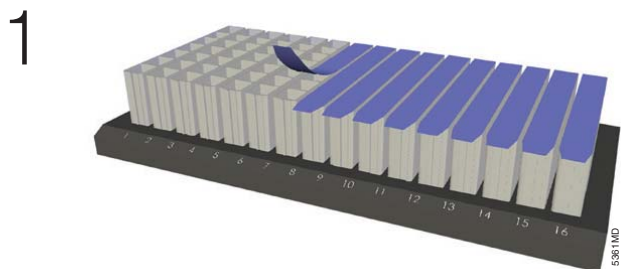




i Please see Maxwell 16 Technical Manual for detailed instructions.

Instructions for Use of Maxwell® 16 Instrument.
Configured in SEV.



Cleaning



Turn off and unplug instrument.



1. Extend platform and wipe off with 70% ethanol then dH₂O.



2. Wipe down magnetic rods and plunger bar adaptor with 70% ethanol.



3. Wipe underside of plunger bar with 70% ethanol then dH₂O.



4. Clean platform slides with a dry cloth.