

Trans-Blot[®] Turbo[™] Instrument



Introduction

This kit provides the necessary tools to help you maintain the pogo pins located in the cavity of your Trans-Blot Turbo instrument. The below procedure should be carried out every six months in order to maintain the performance of your instrument and ensure safety during operation.

Supplied Materials:

- Long cotton swab (15 cm)
- Grease packet

Other Needed Materials:

- Kimwipes
- Isopropyl alcohol
- Laboratory gloves
- Pen or other small blunt object



Maintenance Kit Protocol

Protocol:

Step 1: Turn off and unplug the Trans-Blot Turbo instrument.

Step 2: Wet a kimwipe with isopropyl alcohol and thoroughly wipe each pin in the back of the instrument base cavity. There are eight pins total on each instrument, four per bay.

Step 3: With your gloved finger or the flat end of a pen, push on the end of each pogo pin to ensure they all depress easily. If a pin is stuck, please contact your local Bio-Rad office (details can be found at Bio-Rad.com).

Step 4: Using a second kimwipe, remove excess liquid from each pin.

Step 5: Open the grease packet and place a copious amount on the stick end of the cotton swab.

Step 6: Apply the grease liberally to each pin.

The gold-plated pogo pin electrical contacts on your Trans-Blot Turbo system are now protected from liquid damage and the system is ready for use!

Kimwipes is a trademark of the Kimberly-Clark Corporation.