

Special Instructions to replace anode/cathode electrical contacts on the iBlot® device

Publication Part No. 25-1074

MAN0006751

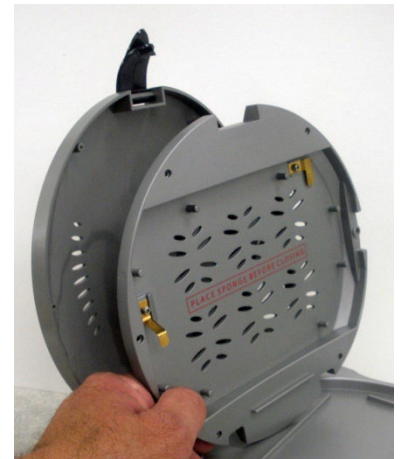
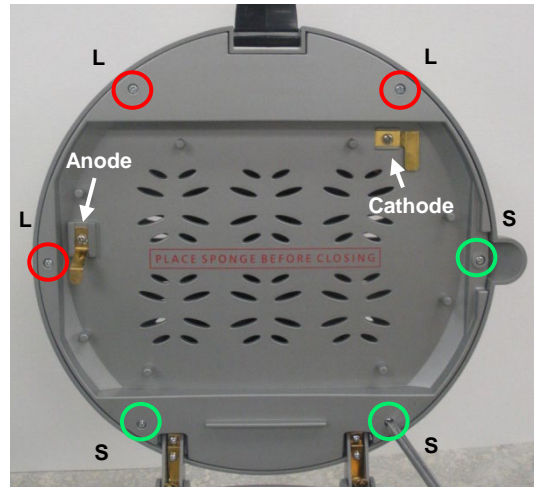
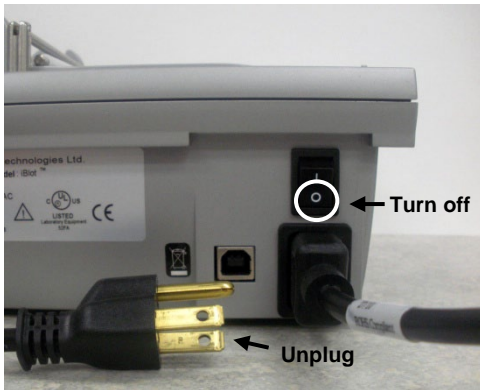
Rev. Date 25 March 2012

Materials required

- Philips-head screwdriver
- Needle-nose pliers
- New iBlot® Device Electrodes (Cat no. IB1002)

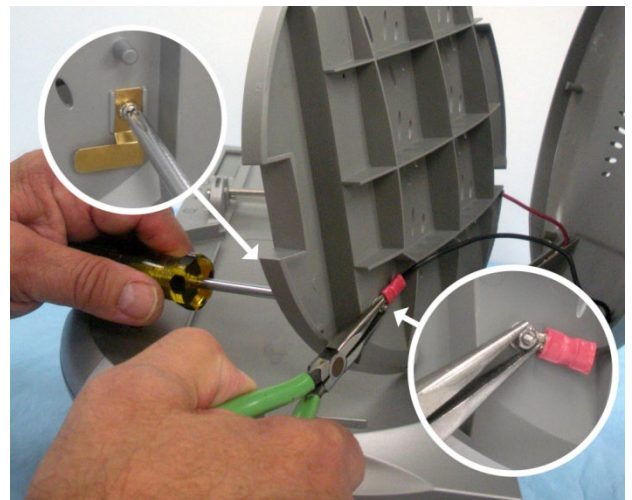
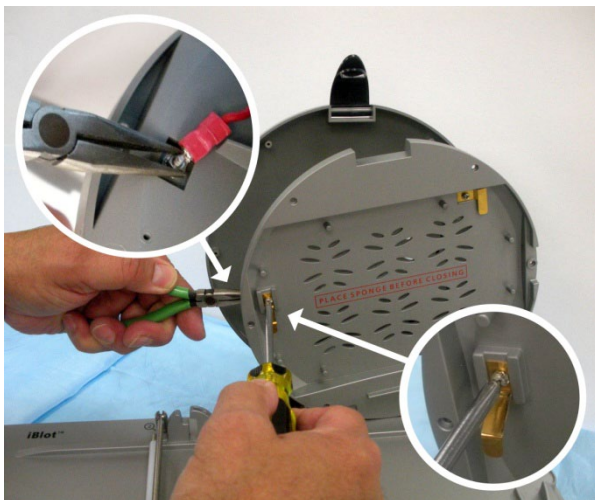
WARNING! Make sure that the iBlot® Gel Transfer Device is turned off and unplugged from any electrical source before replacing the electrodes.

Disassemble lid of iBlot® device



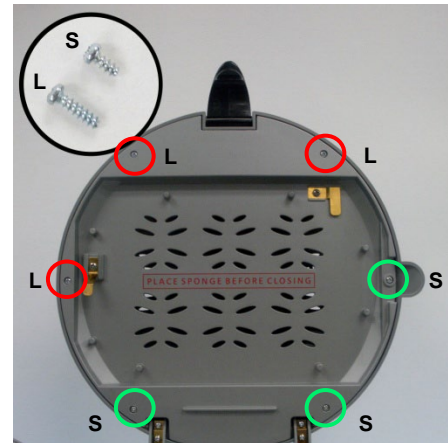
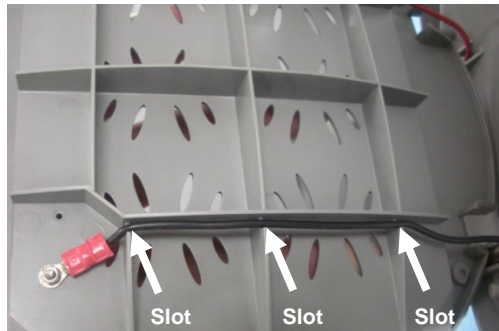
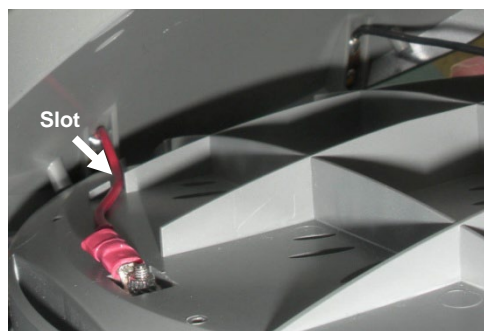
1. Turn off the iBlot® device and unplug the power cord from any electrical source.
2. Use a Philips-head screwdriver to remove the six screws that hold the two portions of the lid together. Place the screws in a secure place. **Note:** There are three long screws (L), and three short screws (S).
3. Separate the two lid components.

Remove electrodes



1. The anode contact is held in place by a nut and screw. Use needle-nose pliers to secure the nut, and then unscrew the screw with the screwdriver to disassemble the contact.
2. The cathode contact is held in place by a nut and screw. Use needle-nose pliers to secure the nut, and then unscrew the screw with the screwdriver to disassemble the contact.

Replace electrodes and reassemble lid of iBlot® device



1. Replace the anode contact, and secure the contact to the wire in reverse order of removal with the appropriate nut and screw.
2. Make sure that the anode wire is tucked into the plastic slot.
3. Replace the cathode contact, and secure the contact to the wire in reverse order of removal with the appropriate nut and screw.
4. Make sure that the cathode wire is tucked into the plastic slots.
5. Reassemble the lid and secure the two lid components with the screws. Remember that the three longer screws (L) are for the top and left holes.

Electrical Safety

Operation of the iBlot® Gel Transfer Device is subject to the following conditions:

- Indoor use
- Altitude below 2000 meters
- Temperature range: 5 to 40°C
- Maximum relative humidity: 80%
- Installation categories (over voltage categories) II; Pollution degree 2
- Mains supply voltage fluctuations not to exceed 10% of the nominal voltage (100–240 V, 50/60 Hz, 3.3 A)
- Mains plug is a disconnect device and must be easily accessible
- Do not attempt to open the iBlot® Gel Transfer Device. To honor the warranty, iBlot® device can only be opened and serviced by Life Technologies
- The protection provided by the equipment may be impaired if the equipment is used in a manner not specified by Life Technologies
- The device must be connected to a mains socket outlet with protective earthing connections
- Ventilation requirements: room ventilation
- The iBlot® Gel Transfer Device complies with part 15 of the FCC rules. Operation of the device is subject to the following two conditions:
 - The device may not cause harmful interference
 - The device must accept any interference received, including interference that may cause undesired operation

Life Technologies Israel Ltd., is the manufacturer and owner of the UL file. For more information, contact:

Life Technologies Israel Ltd.
12 Hamada St.
P.O. Box 4035
Rehovot, Israel 74103

Product Documentation

To obtain a Certificate of Analysis or Safety Data Sheets (SDSs), visit www.lifetechnologies.com/support.

LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

©2012 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their representative owners.

For support visit www.lifetechnologies.com/support or contact techsupport@lifetech.com.

www.lifetechnologies.com

For research use only. Not intended for any animal or human therapeutic or diagnostic use.

Explanation of Certifications, Symbols, and Warnings



Caution



WEEE

The iBlot® Gel Transfer Device complies with the Underwriters Laboratories Inc. regulation, part 15 of the FCC rules, and the European Community Safety requirements. Operation of the iBlot® Gel Transfer Device is subject to the conditions described in this manual. The protection provided by the equipment may be impaired if the equipment is used in a manner not specified by Life Technologies.

The **Caution** symbol denotes a risk of safety hazard. Refer to accompanying documentation to avoid possible personal injury or instrument damage.

The **WEEE** (Waste Electrical and Electronic Equipment) symbol indicates that this product should not be disposed of in unsorted municipal waste. Follow local municipal waste ordinances for proper disposal provisions to reduce the environmental impact of WEEE. Visit <http://www.lifetechnologies.com/weee> for collection and recycling options.

Limited Use Label License: Research Use Only

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, contact outlicensing@lifetech.com or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

Limited Warranty

Life Technologies and/or its affiliate(s) warrant their products as set forth in the Life Technologies General Terms and Conditions of Sale found on the Life Technologies web site at <http://www.lifetechnologies.com/termsandconditions>. If you have any questions, please contact Life Technologies.

life
technologies™