



SERVICE GUIDE

GLH 850 General Laboratory Homogenizer Brush Replacement

Version: E Date: October 2024
Document Number: 03-270BR

Data herein has been verified and validated. It is believed adequate for the intended use of the instrument. If the instrument or procedures are used for purposes over and above the capabilities specified herein, confirmation of the validity and suitability should be obtained; otherwise Revvity does not guarantee results and assumes no obligation or liability. This publication is not a license to operate under, or a recommendation to infringe upon, any process patents.

This product is warranted to be free from defects in material and workmanship for a period of ONE YEAR from the date of delivery. Revvity will repair or replace and return free of charge any part which is returned to its factory within said period, transportation prepaid by user, and which is found upon inspection to have been defective in materials or workmanship. This warranty does not include normal wear from use; it does not apply to any instrument or parts which have been altered by anyone other than an employee of Revvity nor to any instrument which has been damaged through accident, negligence, failure to follow operating instructions, the use of electric currents or circuits other than those specified on the plate affixed to the instrument, misuse, or abuse. Revvity reserves the right to change, alter, modify, or improve any of its instruments without any obligation whatever to make corresponding changes to any instrument previously sold or shipped.

THE FORGOING OBLIGATION IS IN LIEU OF ALL OBLIGATIONS AND LIABILITIES INCLUDING NEGLIGENCE AND ALL WARRANTIES OF MERCHANTABILITY OR OTHERWISE, EXPRESSED OR IMPLIED IN FACT OR BY LAW, AND STATE OUR ENTIRE AND EXCLUSIVE LIABILITY AND BUYERS EXCLUSIVE REMEDY FOR ANY CLAIM OF DAMAGES IN CONNECTION WITH THE SALE OR FURNISHING OF GOODS OR PARTS, THEIR DESIGN, SUITABILITY FOR USE, INSTALLATION, OR OPERATION. REVVITY WILL IN NO EVENT BE LIABLE FOR ANY SPECIAL OR CONSEQUENTIAL DAMAGES WHATSOEVER, AND THEIR LIABILITY UNDER NO CIRCUMSTANCES WILL EXCEED THE CONTRACT PRICE FOR THE GOODS FOR WHICH LIABILITY IS CLAIMED.

This product has been engineered for safety; however, basic safety precautions and common sense must always be demonstrated when using any electrical product.

- Use this product only for its intended purpose.
- Keep this product away from heated surfaces.
- **DO NOT** attempt to modify any part of this product.
- **DO NOT** allow the machine to be submerged in any liquid.
- **DO NOT** use in any setting other than an indoor laboratory.
- **DO NOT** use attachments not recommended by the manufacturer.
- **DO NOT** operate the product if it is damaged in any way.
- **DO NOT** operate the product with the safety ground disconnected.
- **DO NOT** modify the plug or cord that is provided.

Copyright

This manual is published by Revvity 935 Cobb Place Blvd, Kennesaw, GA 30144 USA. Copyright Revvity, Inc. All rights reserved, including but not limited to those rights to reproduce this publication or parts thereof. Reproduction of this publication or parts thereof, or the products it describes, by any means or in any form is expressly prohibited without the written permission of Revvity.

Trademarks

Revvity is a registered trademark of Revvity. All other trademarks are the property of their respective owners.

Procedure

These instructions are written for the replacement of the carbon brushes of the GLH 850 due to wear. Before proceeding ensure the unit is powered off and unplugged.

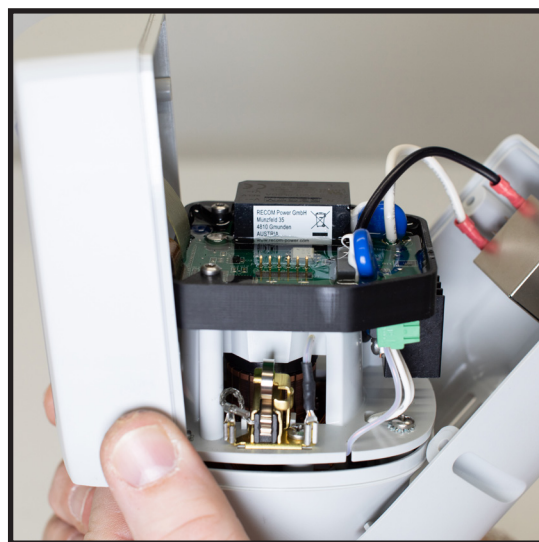
Required equipment and tools.

1. #2 Phillips screwdriver.
2. Small flathead screwdriver.
3. Needle nose pliers.
4. GLH850 Replacement brushes. 110V - 20-110-05R or 220V - 20-220-05R



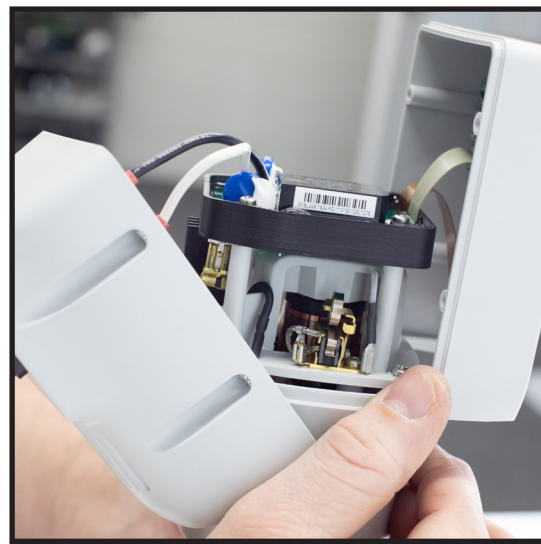
1. Remove the four Phillips screws from the back of the unit.

2. Pull the rear cover straight back. **NOTE:** You do not need to remove the wires from the power block but may do so for convenience. A slight side to side motion while pulling on the connector will aid in removal. Do not pull on the wire.



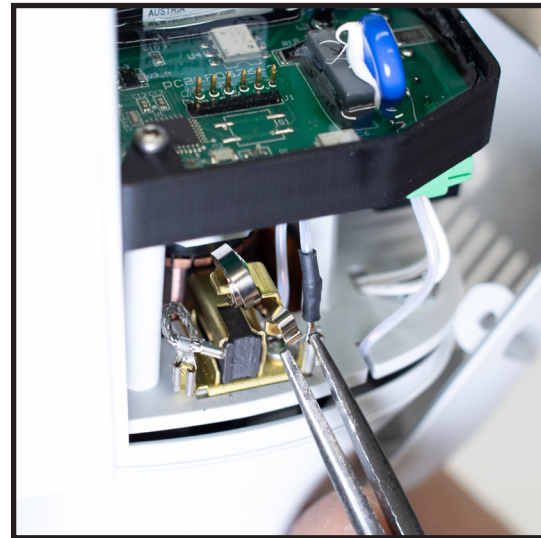
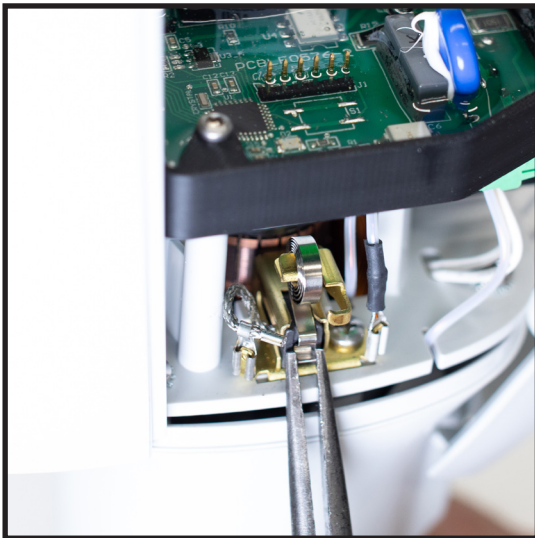
3. Carefully remove the front face. Grasp the unit with both hands on the lower portion of the GLH850 housing. Your thumbs should be placed on the lower edges of the front face, at or below the lower screw holes.

4. Your index fingers should be on the lower portion of the face plate .

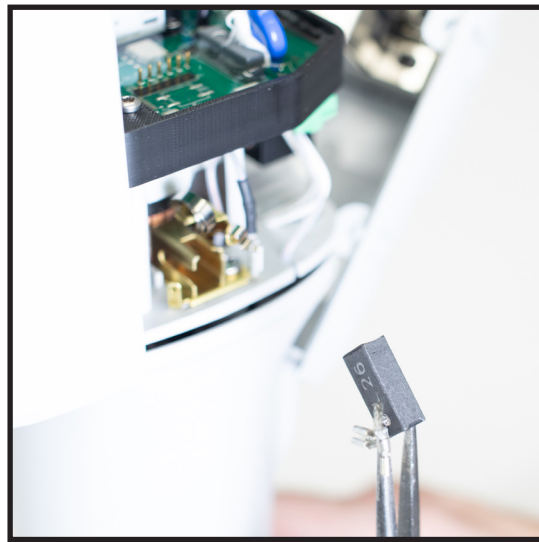
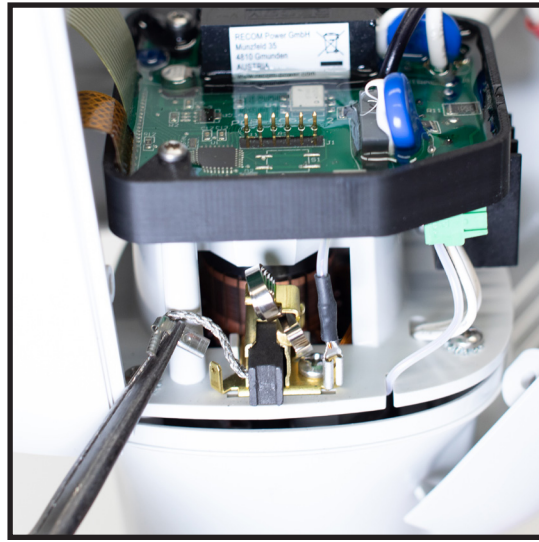
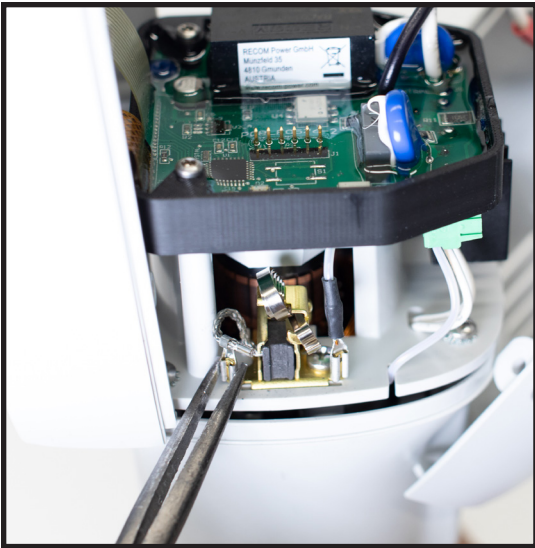


5. Gently push with your thumbs while maintaining slight resistance with your index fingers.

CAUTION: Care must be taken when removing the front face to not damage the control panel and display ribbon cables. These ribbon cables are fragile and can not be repaired. If the ribbon cables are damaged a new face and control board will be required.



6. Using needle nose pliers, carefully place the coil spring onto the coil spring tab of the brush holder.



7. Carefully disconnect the brush lead and remove the brush.

8. Insert new brushes and reset coil spring.

9. Connect brush lead.

10. Re-seat face plate and rear cover.

NOTE: If power wires were disconnected the black wire connects to the top pin and the white wire connects to the bottom.

11. Re-install the four screws.

12. Connect power, turn on unit and verify proper function.

This page was left blank intentionally

This page was left blank intentionally



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.

www.revvity.com

For research use only. Not for use diagnostic procedures

Revvity, Inc.
940 Winter Street
Waltham, MA 02451
USA
www.revvity.com

For a complete listing of our global offices, visit www.revvity.com
Copyright ©2024, Revvity, Inc. All rights reserved.