



## SERVICE GUIDE

# Generator Probes

Version: N

Date: October 2024

Document Number: H258G

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.
- **WARNING:** Always submerge probe in liquid before starting. Never run dry.

**CAUTION:** PTFE lower bearings will wear prematurely if not immersed during operation.

**CAUTION:** Do not attempt to autoclave or lubricate stainless steel bearings.

### **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.

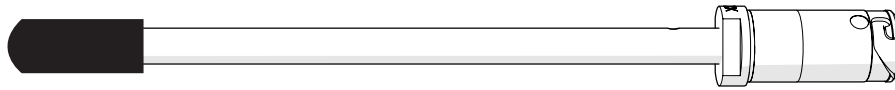
### **Trademarks**

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

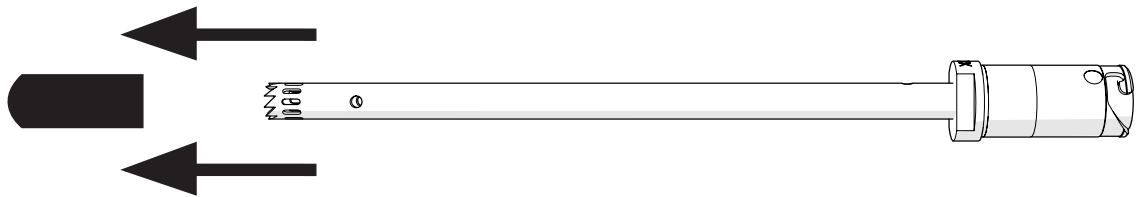
# IMPORTANT INFORMATION

The blue cap must be removed from the end of the generator probe before operating.

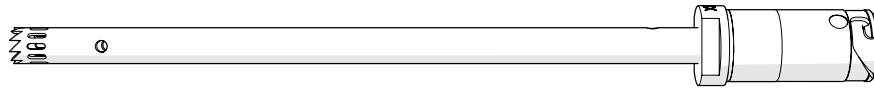
1.



2.



3.

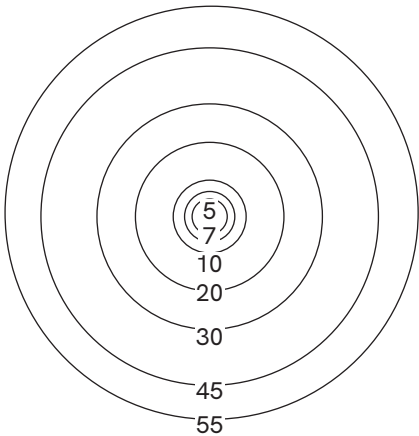


Ready to homogenize.

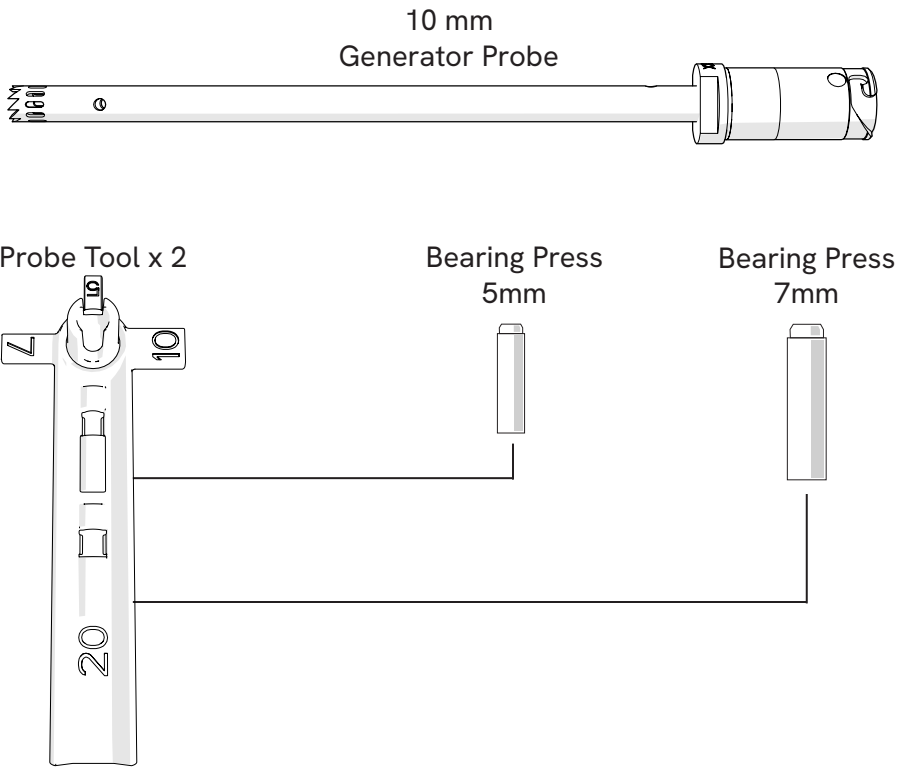
Save the blue cap to protect the probe when it is not in use.

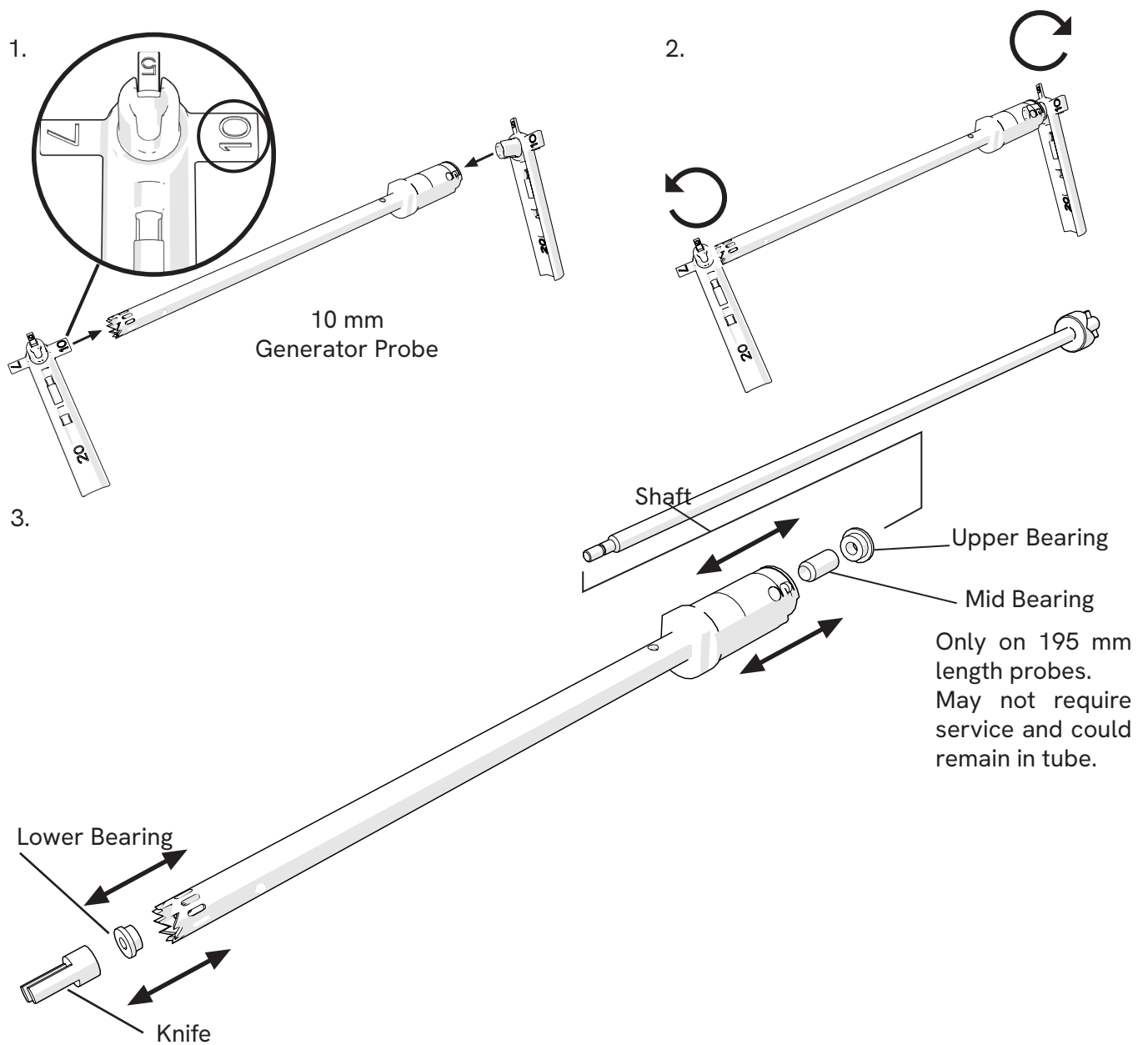
# Generator Probe Size Guide

Match the bottom of the generator probe to determine the diameter in millimeters.



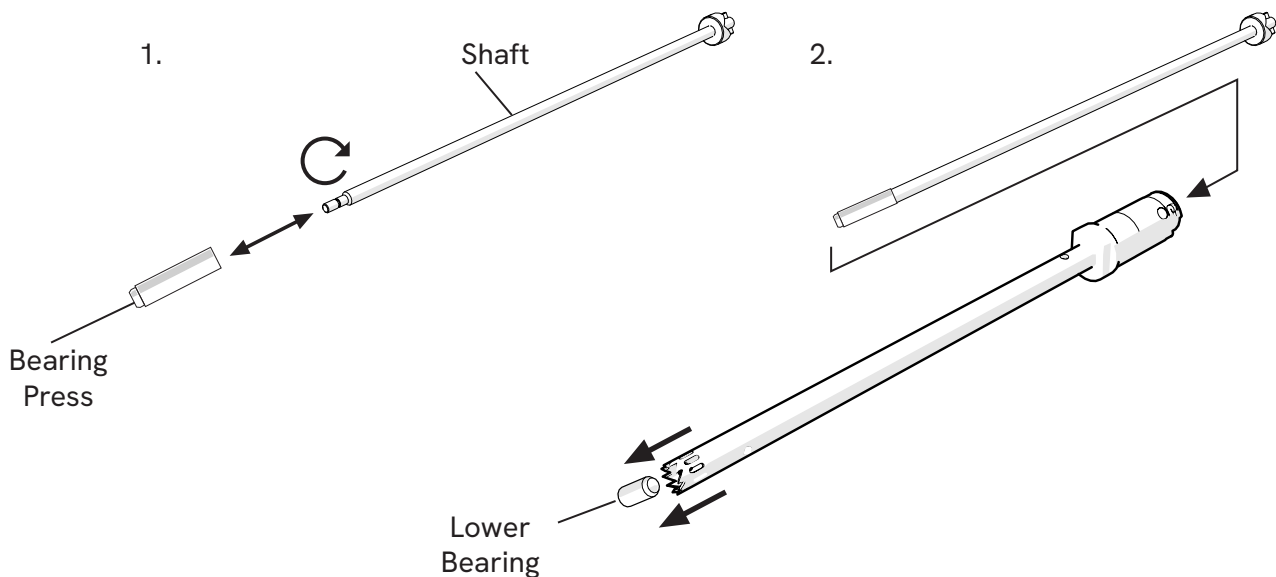
## 5-20 mm Bayonet Style Generator Probes



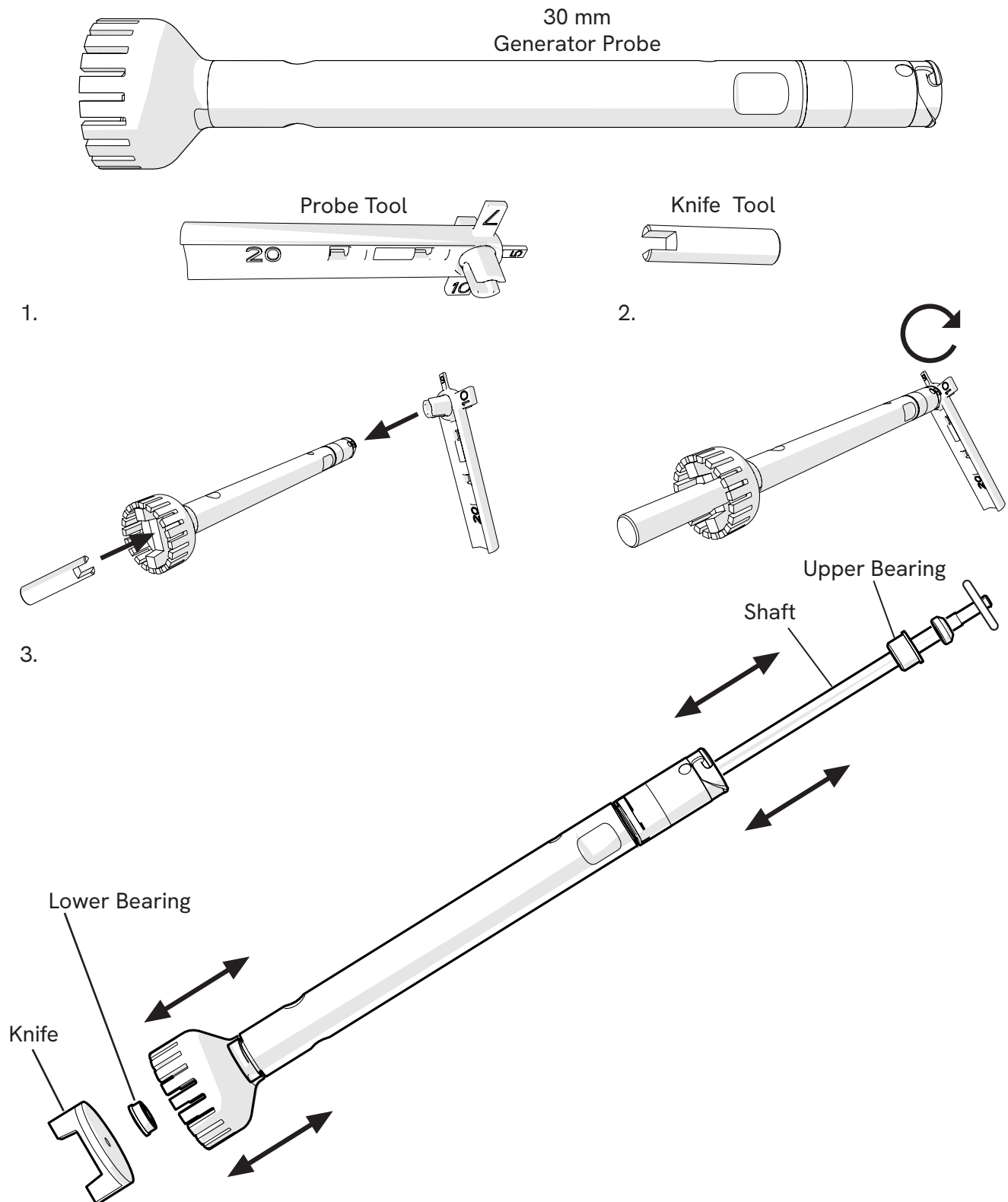


### 5mm and 7mm Generator Probes

1. Attach appropriate Bearing Press to shaft.
2. Insert shaft with attached Bearing Press into probe to remove lower bearing.



## 30-55 mm Bayonet Style Generator Probes



Bearing wear will be determined by factors such as frequency of use, cycle time and processing speed. Long cycle times at high speeds may accelerate bearing wear and require more frequent bearing replacement.

If extended cycle times and/or high speeds are used to homogenize the sample, we recommend changing the upper PTFE bearing to the optional stainless-steel ball bearing. Failure to maintain the bearings may cause damage to your probe.

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](http://www.revvity.com)**

**For research use only. Not for use diagnostic procedures**

Revvity, Inc.  
940 Winter Street  
Waltham, MA 02451  
USA  
[www.revvity.com](http://www.revvity.com)

For a complete listing of our global offices, visit [www.revvity.com](http://www.revvity.com)  
Copyright ©2024, Revvity, Inc. All rights reserved.