

VesselVue™ Microbubble Reagent

P-formulation

Product No.: P-4007003

DESCRIPTION

VesselVue™ formulations are microbubble contrast reagents for contrast enhanced ultrasound (CEUS) imaging and preclinical therapy applications. VesselVue microbubbles can enable the visualization and quantification of tumor or organ vascularity in CEUS imaging, revealing dynamic changes in blood vessels for disease characterization and therapeutic response studies.

The P-formulation (polydisperse) is comprised of microbubbles with a wider size distribution and a smaller mean diameter, making the P-formulation an excellent all-around vascular imaging contrast agent for general CEUS imaging applications.

CONTENTS

- 2 vials – VesselVue Microbubble Contrast Agent, P-formulation
- 1 mL per vial

STORAGE & HANDLING

- Upon receipt, VesselVue P-formulation agents should be IMMEDIATELY STORED AT 4-8°C.
- Do not freeze.
- When stored and handled properly, an unopened (sealed) vial is stable for a minimum of 6 months from date of shipping.
- Allow the probe to warm up to room temperature and activate microbubbles before injecting into an animal.

PROTOCOL

- VesselVue P-formulation must be activated immediately before use with the Bead Ruptor VesselVue Homogenizer, available from Revvity.
- Please scan the QR code on the product packaging or see VesselVue P-formulation User Manual available on our website.

Property	Specification
Shell Type	Phospholipid
Core Type	Perfluorocarbon
Mean Diameter	0.99+/-0.56 µm
Mode Diameter	0.6 µm
Median Diameter	0.83 µm
Size Profile	Poly Distribution
Concentration	1.00 +/- 0.20 x 10 ¹⁰ bubbles/mL
Volume per vial	1 mL
Vial size	3 ml
Shipping condition	Shipped on blue ice

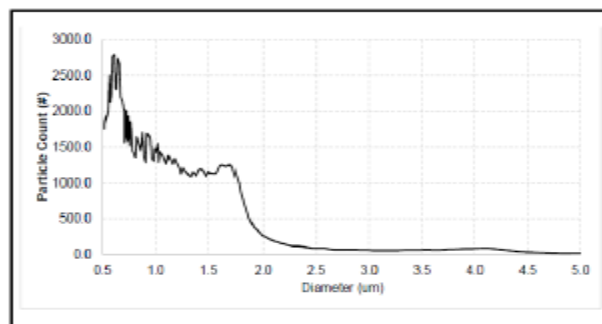


Figure 1: A histogram showing the size distribution of VesselVue P-formulation.

Limited Warranty: This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. 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.

revvity

Revvity
940 Winter Street
Waltham, MA 02451 USA

(800) 762-4000
www.revvity.com

For a complete listing of our global offices, visit www.revvity.com
Copyright ©2023, Revvity, Inc. All rights reserved.