

VesselVue[®] Microbubble Reagent

P-formulation

Product Number: **P-4007001**

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 compared to the S-formulation, 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 VesselVue Microbubble Mixer, available from PerkinElmer.
- Please see In Vivo Imaging Protocol for VesselVue, P-formulation 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
Shipping condition	Shipped on blue ice

