

Product Number:	103526
Lot Number:	403316
Expiration:	2-Feb-2026
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.23×10^9 Beads/mL

Review/Approval

Greg Simone

Quality Events and Training Specialist

[Signature] 09 Oct 2025

Name

Title

Signature/Date