

Product Number:	103527
Lot Number:	309311
Expiration:	3-Apr-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.45 x 10 ⁹ Beads/mL

Review/Approval

Cody Carlson
Name

QC Manager
Title

[Signature] 13 FEB 2024
Signature/Date