

Product Number:	103208
Lot Number:	300911
Expiration:	9-Jan-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.27 x 10 ⁹ Beads/mL

Review/Approval

Brian Ciccarciello
Name

VP, QA+RA
Title

B-C 19 APR 2023
Signature/Date