

Product Number:	103208
Lot Number:	317113
Control Number:	317113A
Expiration:	20-Jun-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.28×10^9 Beads/mL

Review/Approval

Sonia Khillan
Name

Senior QC Analyst
Title

Sonia Khillan 02 NOV 2023
Signature/Date