

Product Number:	103208
Lot Number:	331511
Expiration:	11-Nov-2024
Storage:	2-8°C

Laboratory Analysis		
Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.16 x 10 ⁹ Beads/mL

Review/Approval

Heidi Caswell
Name

Director, QC
Title

Heidi Caswell 01May2024
Signature/Date