

Product Number:	103208
Lot Number:	333513
Expiration:	1-Dec-2024
Storage:	2-8°C

Laboratory Analysis		
Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.29 x 10 ⁹ Beads/mL

Review/Approval

ROBERT SOMOGYI
Name

QC MANAGER
Title

[Signature] 02 MAY 2024
Signature/Date