

Product Number:	103208
Lot Number:	429122
Expiration:	17-Oct-2025
Storage:	2-8°C

Laboratory Analysis		
Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.26 x 10 ⁹ Beads/mL

Review/Approval

Robert Somogyi
Name

QC Manager
Title


Signature/Date

14 NOV 2024