

Product Number:	103208
Lot Number:	429123
Expiration:	17-Oct-2025
Storage:	2-8°C

Laboratory Analysis		
Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.15 x 10 ⁹ Beads/mL

Review/Approval

Robert Somogyi

QC Manager

 05 DEC 2024

Name

Title

Signature/Date