

Product Number:	103208
Lot Number:	429124
Expiration:	17-Oct-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.21 x 10 ⁹ Beads/mL

Review/Approval

Robert Somogyi
Name

QC Manager
Title

 04 Feb 2025
Signature/Date