

Product Number:	103528
Lot Number:	523315
Expiration:	21-Aug-2027
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.32×10^9 Beads/mL

Review/Approval

Kathleen Cornell
Name

Quality Control
Supervisor
Title

[Signature]
04 Dec 2025
Signature/Date