

Product Number:	103526
Lot Number:	302009
Expiration:	20-Jan-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.35×10^9 Beads/mL

Review/Approval

Matt Kaulouris
Name

QC Supervisor
Title

[Signature] 07Feb2023
Signature/Date