

Product Number:	103526
Lot Number:	123211
Expiration:	20-Aug-2022
Storage:	2 - 8°C

Laboratory Analysis

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

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 06Oct2024
Signature/Date