

Product Number:	103526
Lot Number:	202105
Expiration:	21-Jan-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.244 x 10 ⁹ Beads/mL

Review/Approval

Shivani Goel
Name

QC, Manager
Title

Shivani Goel 10 Feb 2022
Signature/Date