

Product Number:	103208
Lot Number:	025510
Expiration:	11-Sep-2021
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.38x10 ⁹ Beads/mL

Review/Approval

Shivani Goel
~~Sivasankari Murugan~~
Name

QC Manager
~~QC Supervisor~~
Title

Shivani Goel 25 Sep 2020
Signature/Date