

Product Number:	103529
Lot Number:	111222
Expiration:	22-Apr-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.49x10 ⁹ Beads/mL

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 03 Sep 2021
Signature/Date