

Product Number:	103526
Lot Number:	129812
Expiration:	25-Oct-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.228 x 10 ⁹ Beads/mL

Review/Approval

Shivani Goel

Name

QC

Title

Shivani Goel 13 Dec 2021

Signature/Date