

Product Number:	103208
Lot Number:	329317
Expiration:	20-Oct-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.27 x 10 ⁹ Beads/mL

Review/Approval

Pratik Shetty
Name

QC
SUPERVISOR
Title

Pratik Shetty
Signature/Date

08 DEC 2023