




Product Number:	103208
Lot Number:	306011
Expiration:	1-Mar-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.31×10^9 Beads/mL

Review/Approval

Name: Prakash Smet Title: QC Supervisor Signature/Date: Pratik Shetty 22 MAY 2023