

Product Number:	103529
Lot Number:	201416
Expiration:	14-Jan-2023
Storage:	2 - 8°C


Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.10 X 10 ⁹ Beads/mL

Review/Approval

Richa Shrestha

QC Supervisor

 21Apr2022

Name

Title

Signature/Date