

Product Number:	103529
Lot Number:	026519
Expiration:	21-Sep-2021
Storage:	2 - 8°C


Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.25x10 ⁹ Beads/mL

Review/Approval

Junpil Park

QC Supervisor

 14 Jan 2021

Name

Title

Signature/Date