

Product Number:	103529
Lot Number:	026517
Expiration:	21-Sep-2021
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.15x10 ⁹ Beads/mL

Review/Approval

Shivani Goel

QC Manager

Shivani Goel ~~SG~~ ^{02 NOV 2020}
Signature/Date 02 NOV 2020

Name

Title

Signature/Date