

Product Number:	103529
Lot Number:	211812
Expiration:	28-Apr-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.49×10^9 Beads/mL

Review/Approval

Matt Smart
Name

QC
Title

MAT 13 Jun 2022
Signature/Date