

Product Number:	103529
Lot Number:	211811
Expiration:	28-Apr-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.33×10^9 Beads/mL

Review/Approval

Matt Smart
Name

Director, QC
Title

MAS 19 May 2022
Signature/Date