

Product Number:	103208
Lot Number:	211206
Expiration:	22-Apr-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.32 x 10 ⁹ Beads/mL

Review/Approval

Matt Smart
Name

QC

Title

[Signature] 17JUN2022
Signature/Date