

Product Number:	103528
Lot Number:	211912
Expiration:	29-Apr-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.59 x 10 ⁹ Beads/mL

Review/Approval

Matt Smart
Name

QC
Title

M SA 025m 2022
Signature/Date