

Product Number:	103208
Lot Number:	221510
Expiration:	3-Aug-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.21 x 10 ⁹ Beads/mL

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan / 04 OCT 2022
Signature/Date