

Product Number:	103208
Lot Number:	230807
Expiration:	4-Nov-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.26 x 10 ⁹ Beads/mL

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan / 11 JAN 2023
Signature/Date