

Product Number:	103527
Lot Number:	222008
Expiration:	8-Aug-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.34 x 10 ⁹ Beads/mL

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan / 22 DEC 2022
Signature/Date