

Product Number:	103528
Lot Number:	211914
Expiration:	29-Apr-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.18 x 10 ⁹ Beads/mL

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan / 07NOV2022
Signature/Date