

Product Number:	103528
Lot Number:	303310
Expiration:	2-Feb-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.32 x 10 ⁹ Beads/mL

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan 25 MAY 2023
Signature/Date