

Product Number:	103529
Lot Number:	231107
Expiration:	7-Nov-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.28 x 10 ⁹ Beads/mL

Review/Approval

Matt Koudouris
Name

QC Supervisor
Title

[Signature]
Signature/Date