

Product Number:	103529
Lot Number:	226222
Expiration:	19-Sep-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.40 x 10 ⁹ Beads/mL

Review/Approval

Matt Koubouris

QC Supervisor

[Signature] 24 Jan 2023

Name

Title

Signature/Date