

Product Number:	103208
Lot Number:	231918
Expiration:	15-Nov-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.17 x 10 ⁹ Beads/mL

Review/Approval

Sonia Killian

Name

QC Manager

Title

Sonia Killian 09MAR2023

Signature/Date