

Product Number:	103208
Lot Number:	331509
Expiration:	11-Nov-2024
Storage:	2-8°C

Laboratory Analysis

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

Review/Approval

ROBERT SOMOGYI
Name

QC MANAGER
Title

[Signature] 19 JAN 2024
Signature/Date