

Product Number:	103208
Lot Number:	334715
Expiration:	13-Dec-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.16 x 10 ⁹ Beads/mL


Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title

 27 FEB 2024

Signature/Date