

Product Number:	103208
Lot Number:	406007
Expiration:	28-Feb-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.23 x 10 ⁹ Beads/mL


Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title

 25 MAR 2024

Signature/Date