

Product Number:	103208
Lot Number:	407913
Expiration:	19-Mar-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.26 x 10 ⁹ Beads/mL

Review/Approval

Robert Somogyi

QC Manager

 05 JUN 2024

Name

Title

Signature/Date