

Product Number:	103208
Lot Number:	411740
Expiration:	26-Apr-2025
Storage:	2-8°C

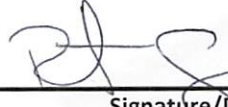
Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.25 x 10 ⁹ Beads/mL

Review/Approval

Robert Somogyi
Name

QC Manager
Title

 24 JUL 2024
Signature/Date