

Product Number:	103208
Lot Number:	505814
Expiration:	27-Feb-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.27×10^9 Beads/mL

Review/Approval

Robert Somogyi

Name

QC Manager

Title

 22 APR 2015
Signature/Date