

Product Number:	103612
Lot Number:	209716
Expiration:	5-Dec-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	3.26×10^9 Beads/mL

Review/Approval

Matt Smart

Name

QC

Title

MAA 10 May 2022

Signature/Date