

Product Number:	103527
Lot Number:	210513
Expiration:	15-Apr-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.45×10^9 Beads/mL

Review/Approval

Matt Smart

Name

QC

Title

MAT 02 Jun 2022

Signature/Date