

Product Number:	103526
Lot Number:	210514
Expiration:	15-Apr-2023
Storage:	2 - 8°C

Laboratory Analysis

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

Review/Approval

Matt Smart

Name

QC

Title

Matt Smart 30 Jun 2022

Signature/Date