

Product Number:	103529
Lot Number:	213211
Expiration:	12-May-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.18 x 10 ⁹ Beads/mL

Review/Approval

Matt Smart

Name

QC

Title

MA 26 Sep 2022

Signature/Date