

Product Number:	100993
Lot Number:	222809
Expiration:	16-Aug-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	5.20 x 10 ⁶ Beads/mL

Review/Approval

Matt Smart
Name

QC
Title

MSA 22Aug2022
Signature/Date