

Product Number:	100998
Lot Number:	211109
Expiration:	21-Apr-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	PASS

Total Bead Concentration (beads/mL)	5.41 x 10 ⁶
--	------------------------

Review/Approval

Richa Shrestha
Name

QC Supervisor
Title

Richa Shrestha 29 Apr 2023
Signature/Date