

Product Number:	103528
Lot Number:	100826
Expiration:	8-Jan-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.17×10^9 Beads/mL

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 06 May 2021
Signature/Date