

Product Number:	103526
Lot Number:	304415
Expiration:	13-Feb-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.27×10^9 Beads/mL

Review/Approval

Sonia Khilken

Name

QC Manager

Title

Sonia Khilken 13 OCT 2023

Signature/Date