

Product Number:	103526
Lot Number:	007212
Expiration:	12-Mar-2021
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.36x10 ⁹ Beads/mL

Review/Approval

Patrizia Stadler

Name

QC Assoc. Director

Title

P. Stadler

Signature/Date

01 May 2020