

Product Number:	104006
Lot Number:	328215
Expiration:	9-Oct-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.31 x 10 ⁹ Beads/mL

Review/Approval

Sonia Khillan	Senior QC Analyst	Sonia Khillan 25 Oct 2023
Name	Title	Signature/Date