

Product Number:	103612
Lot Number:	209719
Expiration:	5-Dec-2022
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	3.23 x 10 <sup>9</sup> Beads/mL

Review/Approval

*Matt Smart*

Name

QC

Title

*MRT 27 Jun 2022*

Signature/Date