

Product Number:	101076
Lot Number:	503106
Expiration:	13-Jul-2022

Component	Part Number	Lot Number	Bead Count (beads/mL)
Carboxylated Paramagnetic Beads	103611	102534	3.05 x 10 ⁹

Component	Part Number	Lot Number	Concentration (nM)
SBG Concentrate	103397	119416	50

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 17 Aug 2021
Signature/Date

Sta _____ AS.