

Product Number:	101076
Lot Number:	503250
Expiration:	13-Jul-2022

Component	Part Number	Lot Number	Bead Count (beads/mL)
Carboxylated Paramagnetic Beads	103611	108822	2.49 x 10 ⁹

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

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 07Dec2021
Signature/Date

Statement: The above product was tested using Quanterix procedures and conforms to Quanterix's specifications.