

Product Number:	101076
Lot Number:	501298
Expiration:	17-Apr-2019

Component	Part Number	Lot Number	Bead Count (beads/mL)
Carboxylated Paramagenetic Beads	103206	MF-052418-01	3.23 x 10 ⁹

Component	Part Number	Lot Number	Concentration (nM)
SBG Concentrate	100439	81072	186

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

 22 Jun 2018
Signature/Date

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