

Product Number:	101076
Lot Number:	501600
Expiration:	29-Aug-2019

Component	Part Number	Lot Number	Bead Count (beads/mL)
Carboxylated Paramagenetic Beads	103206	823619	3.29x10 ⁹

Component	Part Number	Lot Number	Concentration (nM)
SBG Concentrate	100439	824115	155

Review/Approval

Patrizia Stadler

Associate Director,
QC

P. Stadler 23 Jan 2019

Name

Title

Signature/Date

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