

Product Number:	101076
Lot Number:	502122
Expiration:	5-Aug-2020

Component	Item Number	Lot Number	Bead Count (beads/mL)
<b>Carboxylated Paramagenetic Beads</b>	103206	823620	3.41 x 10 <sup>9</sup>

Component	Part Number	Lot Number	Concentration (nM)
<b>SBG Concentrate</b>	103397	921713	50

Review/Approval

Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	19 Nov 2019
Name	Title	Signature/Date	

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