

Product Number:	101076
Lot Number:	502180
Expiration:	5-Aug-2020

Component	Part Number	Lot Number	Bead Count (beads/mL)
Carboxylated Paramagenetic Beads	103206	823622	3.62 x 10 <sup>9</sup>

Component	Part Number	Lot Number	Concentration (nM)
SBG Concentrate	103397	921713	50

Review/Approval Patrizia Stadler Assoc. Director, QC *p-stadler* 30 Dec 2019  
 Name Title Signature/Date

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