

<b>Product Number:</b>	101354
<b>Lot Number:</b>	501670
<b>Expiration:</b>	5-Nov-2019

Component	Item Number	Lot Number
Beads	103207	822906
Detector/Sample Diluent	101359	902927
Biotinylation Reaction Buffer	101358	904302
Bead Conjugation Buffer	101357	830909
Bead Blocking Buffer	101356	830905
Bead Wash Buffer	101355	830904
Bead Diluent Buffer	100458	833308

Bead Count (beads/mL)	3.04x10 <sup>9</sup>
--------------------------	----------------------

Review/Approval

Patrizia Stadler	QC Assoc. Director	<i>P. Stodlev 21Feb2019</i>
Name	Title	Signature/Date

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