

Product Number:	101354
Lot Number:	501780
Expiration:	5-Nov-2019

Component	Item Number	Lot Number
Beads	103207	822906
Detector/Sample Diluent	101359	905819
Biotinylation Reaction Buffer	101358	904308
Bead Conjugation Buffer	101357	830909
Bead Blocking Buffer	101356	906502
Bead Wash Buffer	101355	903616
Bead Diluent Buffer	100458	904217

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

Review/Approval

Patrizia Stadler QC Assoc. Director

*P. Stadler* 22 Apr 2019

Name

Title

Signature/Date

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