

Product Number:	101354
Lot Number:	501852
Expiration:	11-Feb-2020

Component	Item Number	Lot Number
Beads	103207	912111
Detector/Sample Diluent	101359	913112
Biotinylation Reaction Buffer	101358	904308
Bead Conjugation Buffer	101357	916503
Bead Blocking Buffer	101356	906502
Bead Wash Buffer	101355	910013
Bead Diluent Buffer	100458	904223

Bead Count (beads/mL)	3.20x10 <sup>9</sup>
--------------------------	----------------------

Review/Approval

Patrizia Stadler QC Assoc. Director

Name

Title

Signature/Date

*P. Stadler* 26JUN2019

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