

Product Number:	101354
Lot Number:	501473
Expiration:	14-May-2019


Component	Item Number	Lot Number
Beads	103207	822906
Detector/Sample Diluent	101359	827612
Biotinylation Reaction Buffer	101358	MF-051718-07
Bead Conjugation Buffer	101357	MF-051718-05
Bead Blocking Buffer	101356	MF-051418-08
Bead Wash Buffer	101355	MF-051418-10
Bead Diluent Buffer	100458	MF-052318-03

Bead Count (beads/mL)	3.04 x 10 ⁹
--------------------------	------------------------

Review/Approval

Sivasankari Murugan

QC Supervisor

 25 OCT 2018

Name

Title

Signature/Date

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