

Product Number:	101354
Lot Number:	501111
Expiration:	16-Aug-2018

Component	Item Number	Lot Number
Bead Diluent Buffer	100458	MF-081617-03
Bead Wash Buffer	101355	MF-081817-11
Bead Blocking Buffer	101356	MF-081617-09
Bead Conjugation Buffer	101357	MF-081617-06
Biotinylation Reaction Buffer	101358	MF-081817-06
Detector/Sample Diluent	101359	MF-081817-09
Beads	101360	MF-012418-08

Bead Count (beads/mL)	2.80 x 10 ⁹
--------------------------	------------------------

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

 06 FEB 2018
Signature/Date

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