

Product Number:	101354
Lot Number:	501165
Expiration:	16-Aug-2018

Component	Item Number	Lot Number
Beads	101360	MF-030218-03
Detector/Sample Diluent	101359	MF-112417-11
Biotinylation Reaction Buffer	101358	MF-112817-13
Bead Conjugation Buffer	101357	MF-112417-05
Bead Blocking Buffer	101356	MF-112417-07
Bead Wash Buffer	101355	MF-112417-10
Bead Diluent Buffer	100458	MF-081617-03

Bead Count (beads/mL)	2.64 x 10 ⁹
--------------------------	------------------------

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title


Signature/Date 09 MAR 2018

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