

Product Number:	101354
Lot Number:	501282
Expiration:	24-Nov-2018

Component	Item Number	Lot Number
Beads	103207	MF-050718-07
Detector/Sample Diluent	101359	MF-042318-06
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-112817-23

Bead Count (beads/mL)	3.28 x 10 ⁹
--------------------------	------------------------

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title


Signature/Date 12 Jun 2018

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