

Product Number:	101354
Lot Number:	501499
Expiration:	17-May-2019

Component	Item Number	Lot Number
Beads	103207	822906
Detector/Sample Diluent	101359	831007
Biotinylation Reaction Buffer	101358	MF-051718-07
Bead Conjugation Buffer	101357	MF-051718-05
Bead Blocking Buffer	101356	830905
Bead Wash Buffer	101355	830904
Bead Diluent Buffer	100458	830608

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

Review/Approval

Sivasankari Murugan

QC Supervisor

Name

Title

Signature/Date



15 Nov 2018

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