

Product Number:	102040
Lot Number:	501436
Expiration:	14-May-2019

Component	Item Number	Lot Number
Detector/Sample Diluent	101359	823617
Bead Conjugation Buffer	101357	MF-051718-05
Bead Diluent Buffer	100458	MF-052318-03
Biotinylation Reaction Buffer	101358	MF-051718-07
Bead Blocking Buffer	101356	MF-051418-08
Bead Wash Buffer	101355	MF-051418-10

Review/Approval

Sivasankari Murugan QC Supervisor

[Signature] 12 OCT 2018

Name

Title

Signature/Date

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