

Product Number:	102040
Lot Number:	501316
Expiration:	23-Apr-2019

Component	Item Number	Lot Number
Detector/Sample Diluent	101359	MF-042318-06
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  
Name

QC Supervisor  
Title

  
Signature/Date 11/20/2018

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