

Product Number:	102040
Lot Number:	501537
Expiration:	17-May-2019

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

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

  
Signature/Date

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