

Product Number:	102040
Lot Number:	501283
Expiration:	24-Nov-2018

Component	Item Number	Lot Number
Detector/Sample Diluent	101359	MF-042318-06
Bead Conjugation Buffer	101357	MF-112417-05
Bead Diluent Buffer	100458	MF-112817-23
Biotinylation Reaction Buffer	101358	MF-112817-13
Bead Blocking Buffer	101356	MF-112417-07
Bead Wash Buffer	101355	MF-112417-10

Review/Approval

Sivasankari Murugan

QC Supervisor

Name

Title

Signature/Date



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