

Product Number:	102040
Lot Number:	501807
Expiration:	5-Nov-2019

Component	Item Number	Lot Number
Detector/Sample Diluent	101359	905819
Bead Conjugation Buffer	101357	830909
Bead Diluent Buffer	100458	904223
Biotinylation Reaction Buffer	101358	904308
Bead Blocking Buffer	101356	906502
Bead Wash Buffer	101355	910013

Review/Approval

Patrizia Stadler

QC Assoc Director

P. Stadler 03 Nov 2019

Name

Title

Signature/Date

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