

Product Number:	102040
Lot Number:	501652
Expiration:	2-Nov-2019

Component	Item Number	Lot Number
Detector/Sample Diluent	101359	831007
Bead Conjugation Buffer	101357	830909
Bead Diluent Buffer	100458	830608
Biotinylation Reaction Buffer	101358	904302
Bead Blocking Buffer	101356	830905
Bead Wash Buffer	101355	830904

Review/Approval

Patrizia Stadler

QC Assoc
Director

P. Stadler 12 Feb 2019

Name

Title

Signature/Date

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