

Product Number:	101354
Lot Number:	502139
Expiration:	6-Mar-2020

Component	Item Number	Lot Number
Beads	103207	912113
Detector/Sample Diluent	101359	928214
Biotinylation Reaction Buffer	101358	924108
Bead Conjugation Buffer	101357	926906
Bead Blocking Buffer	101356	906502
Bead Wash Buffer	101355	919220
Bead Diluent Buffer	100458	921319

Bead Count (beads/mL)	3.20x10 ⁹
--------------------------	----------------------

Review/Approval

Patrizia Stadler QC Assoc. Director

P. Stadler 26 Nov 2019

Name

Title

Signature/Date

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