

Product Number:	101354
Lot Number:	502654
Expiration:	28-Aug-2021

Component	Item Number	Lot Number
Beads	103612	024813
Detector/Sample Diluent	101359	031411
Biotinylation Reaction Buffer	101358	025909
Bead Conjugation Buffer	101357	034002
Bead Blocking Buffer	101356	024106
Bead Wash Buffer	101355	031113
Bead Diluent Buffer	100458	031128

Bead Count (beads/mL)	3.28x10 ⁹
--------------------------	----------------------

Review/Approval

Junpil Park
Name

QC Supervisor
Title

 14 Dec 2020
Signature/Date

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