

Product Number:	101354
Lot Number:	502677
Expiration:	28-Aug-2021

Component	Item Number	Lot Number
Beads	103612	031523
Detector/Sample Diluent	101359	031906
Biotinylation Reaction Buffer	101358	025909
Bead Conjugation Buffer	101357	034002
Bead Blocking Buffer	101356	024106
Bead Wash Buffer	101355	031113
Bead Diluent Buffer	100458	026620

Bead Count (beads/mL)	3.08x10 ⁹
--------------------------	----------------------

Review/Approval

Junpil Park

QC Supervisor

[Signature] 05 Jan 2021

Name

Title

Signature/Date

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