

Product Number:	101354
Lot Number:	502826
Expiration:	28-Aug-2021

Component	Item Number	Lot Number
Beads	103612	031526
Detector/Sample Diluent	101359	106318
Biotinylation Reaction Buffer	101358	100825
Bead Conjugation Buffer	101357	100703
Bead Blocking Buffer	101356	024111
Bead Wash Buffer	101355	107721
Bead Diluent Buffer	100458	105417

Bead Count (beads/mL)	2.80x10 ⁹
--------------------------	----------------------

Review/Approval

Shivani Goel
Name

QC Manager Shivani Goel 30 Mar 2021
Title Signature/Date

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