

Product Number:	101354
Lot Number:	502881
Expiration:	28-Aug-2021

Component	Item Number	Lot Number
Beads	103612	108824
Detector/Sample Diluent	101359	108821
Biotinylation Reaction Buffer	101358	100825
Bead Conjugation Buffer	101357	100710
Bead Blocking Buffer	101356	024111
Bead Wash Buffer	101355	107721
Bead Diluent Buffer	100458	105417

Bead Count (beads/mL)	3.26x10 ⁹
--------------------------	----------------------

Review/Approval

Junpil Park
Name

QC Supervisor
Title

 17 Apr 2021
Signature/Date

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