

Product Number:	101354
Lot Number:	502631
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	028403
Bead Blocking Buffer	101356	024106
Bead Wash Buffer	101355	031113
Bead Diluent Buffer	100458	031128

Bead Count (beads/mL)	3.28x10 <sup>9</sup>
--------------------------	----------------------

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

*Shivani Goel* 30 NOV 2020  
Signature/Date

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