

Product Number:	101354
Lot Number:	503086
Expiration:	29-Mar-2022

Component	Item Number	Lot Number
Beads	103612	115510
Detector/Sample Diluent	101359	117115
Biotinylation Reaction Buffer	101358	117222
Bead Conjugation Buffer	101357	108832
Bead Blocking Buffer	101356	110523
Bead Wash Buffer	101355	117223
Bead Diluent Buffer	100458	118710

Bead Count (beads/mL)	2.696x10 <sup>9</sup>
--------------------------	-----------------------

Review/Approval

Shivani Goel  
Name

QC  
Title

Shivani Goel  
Signature/Date

03 Aug  
2021

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