

Product Number:	101354
Lot Number:	502900
Expiration:	7-Jan-2022

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	110521
Bead Wash Buffer	101355	107721
Bead Diluent Buffer	100458	105417

Bead Count (beads/mL)	3.26x10 <sup>9</sup>
--------------------------	----------------------

Review/Approval

*Shivani Goel*

Name

*QC Manager*

Title

*Shivani Goel*

Signature/Date

*28 Apr 2021*

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