

Product Number:	101354
Lot Number:	503114
Expiration:	29-Mar-2022

Component	Item Number	Lot Number
Beads	103612	115515
Detector/Sample Diluent	101359	120409
Biotinylation Reaction Buffer	101358	120909
Bead Conjugation Buffer	101357	108832
Bead Blocking Buffer	101356	110523
Bead Wash Buffer	101355	117223
Bead Diluent Buffer	100458	116408

Bead Count (beads/mL)	2.67 x 10 ⁹
--------------------------	------------------------

Review/Approval

Shivani Goel
Name

QC

Title

Shivani Goel 23 Aug 2021
Signature/Date

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