

Product Number:	101354
Lot Number:	502729
Expiration:	28-Aug-2021

Component	Item Number	Lot Number
Beads	103612	031525
Detector/Sample Diluent	101359	102708
Biotinylation Reaction Buffer	101358	100818
Bead Conjugation Buffer	101357	100703
Bead Blocking Buffer	101356	024109
Bead Wash Buffer	101355	031127
Bead Diluent Buffer	100458	035807

Bead Count (beads/mL)	2.99x10 ⁹
--------------------------	----------------------

Review/Approval

Shivani Goel

 Name Title Signature/Date

Shivani Goel 08 Feb 2021

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