


Product Number:	101354
Lot Number:	502689
Expiration:	28-Aug-2021

Component	Item Number	Lot Number
Beads	103612	031523
Detector/Sample Diluent	101359	031906
Biotinylation Reaction Buffer	101358	034007
Bead Conjugation Buffer	101357	034002
Bead Blocking Buffer	101356	024109
Bead Wash Buffer	101355	031127
Bead Diluent Buffer	100458	035807

Bead Count (beads/mL)	3.08x10 ⁹
--------------------------	----------------------

Review/Approval

Junpil Park QC Supervisor  13Jan 2021

 Name Title Signature/Date

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