

Product Number:	101354
Lot Number:	502597
Expiration:	24-May-2021

Component	Item Number	Lot Number
Beads	103612	015313
Detector/Sample Diluent	101359	028407
Biotinylation Reaction Buffer	101358	023711
Bead Conjugation Buffer	101357	028403
Bead Blocking Buffer	101356	024106
Bead Wash Buffer	101355	026618
Bead Diluent Buffer	100458	014512

Bead Count (beads/mL)	3.24x10 ⁹
--------------------------	----------------------

Review/Approval

Shivani Goel

QC Manager

Shivani Goel
Signature/Date

30 Oct 2020

Name

Title

Signature/Date

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