

<b>Product Number:</b>	101354
<b>Lot Number:</b>	502585
<b>Expiration:</b>	26-Apr-2021

Component	Item Number	Lot Number
Beads	103612	015313
Detector/Sample Diluent	101359	028407
Biotinylation Reaction Buffer	101358	025909
Bead Conjugation Buffer	101357	011709
Bead Blocking Buffer	101356	024106
Bead Wash Buffer	101355	026618
Bead Diluent Buffer	100458	022412

Bead Count (beads/mL)	3.24x10 <sup>9</sup>
--------------------------	----------------------

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

*Shivani Goel 2004 2020*  
Signature/Date

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