

Product Number:	101354
Lot Number:	503438
Expiration:	2-Oct-2022

Component	Item Number	Lot Number
488-dyed Singleplex Beads <sup>1</sup>	104006	205909
Detector/Sample Diluent	101359	206616
Biotinylation Reaction Buffer	101358	209003
Bead Conjugation Buffer	101357	134611
Bead Blocking Buffer	101356	127515
Bead Wash Buffer	101355	133220
Bead Diluent Buffer	100458	135012

<sup>1</sup> Assay Beads provided are conjugated with 488 nm fluorescent dye  
For more information, please visit

<https://go.quanterix.com/HomebrewBeadUpdate>



Bead Count (beads/mL)	1.17 x 10 <sup>9</sup>
--------------------------	------------------------

Review/Approval

Shivani Goel

QC Manager

*Shivani Goel* 11Apr2022

Name

Title

Signature/Date

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