

Product Number:	101354
Lot Number:	503348
Expiration:	27-Jul-2022

Component	Item Number	Lot Number
488-dyed Singleplex Beads <sup>1</sup>	104006	131609
Detector/Sample Diluent	101359	134925
Biotinylation Reaction Buffer	101358	120910
Bead Conjugation Buffer	101357	134608
Bead Blocking Buffer	101356	127515
Bead Wash Buffer	101355	133218
Bead Diluent Buffer	100458	134609

<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)	1.31 x 10 <sup>9</sup>
--------------------------	------------------------

Review/Approval

Shivani Goel

QC, Manager

*Shivani Goel* 10Feb2022

Name

Title

Signature/Date

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