

Product Number:	101354
Lot Number:	503447
Expiration:	2-Oct-2022

Component	Item Number	Lot Number
488-dyed Singleplex Beads <sup>1</sup>	104006	131517
Detector/Sample Diluent	101359	134921
Biotinylation Reaction Buffer	101358	209003
Bead Conjugation Buffer	101357	134608
Bead Blocking Buffer	101356	127513
Bead Wash Buffer	101355	127517
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)	1.33 x 10 <sup>9</sup>
--------------------------	------------------------

Review/Approval

Richa Shrestha  
Name

QC Supervisor  
Title

*[Signature]*  
Signature/Date  
15AIX2022

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