

Product Number:	101354
Lot Number:	503321
Expiration:	27-Jul-2022

Component	Item Number	Lot Number
488-dyed Singleplex Beads ¹	104006	131517
Detector/Sample Diluent	101359	134921
Biotinylation Reaction Buffer	101358	120910
Bead Conjugation Buffer	101357	134608
Bead Blocking Buffer	101356	127513
Bead Wash Buffer	101355	127517
Bead Diluent Buffer	100458	122710

¹ 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 ⁹
--------------------------	------------------------

Review/Approval

Ruby Burke
Name

QC, Sr. Director
Title

Ruby Burke 25Jan2022
Signature/Date

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