

Product Number:	101354
Lot Number:	503791
Expiration:	6-Oct-2023

Component	Item Number	Lot Number
488-dyed Singleplex Beads ¹	104006	231420
Detector/Sample Diluent	101359	300910
Biotinylation Reaction Buffer	101358	234206
Bead Conjugation Buffer	101357	233907
Bead Blocking Buffer	101356	227910
Bead Wash Buffer	101355	234302
Bead Diluent Buffer	100458	234304

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)	1.22 x 10 ⁹
--------------------------	------------------------

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Soni Kelle 25JAN2023
Signature/Date