

Product Number:	101354
Lot Number:	503690
Expiration:	2-Mar-2023

Component	Item Number	Lot Number
488-dyed Singleplex Beads ¹	104006	222307
Detector/Sample Diluent	101359	218620
Biotinylation Reaction Buffer	101358	220915
Bead Conjugation Buffer	101357	223710
Bead Blocking Buffer	101356	206118
Bead Wash Buffer	101355	206115
Bead Diluent Buffer	100458	223708

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 QC Manager Sonia Khillan / 08SEP2022
Name Title Signature/Date