

Product Number:	101354
Lot Number:	503716
Expiration:	25-Aug-2023

Component	Item Number	Lot Number
488-dyed Singleplex Beads ¹	104006	226507
Detector/Sample Diluent	101359	227312
Biotinylation Reaction Buffer	101358	227604
Bead Conjugation Buffer	101357	227705
Bead Blocking Buffer	101356	227617
Bead Wash Buffer	101355	227908
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)	1.32 x 10 ⁹
--------------------------	------------------------

Review/Approval

Matt Koulouris
Name

QC Supervisor
Title

[Signature] 11 Oct 2022
Signature/Date