

Product Number:	101354
Lot Number:	503506
Expiration:	2-Mar-2023

Component	Item Number	Lot Number
488-dyed Singleplex Beads ¹	104006	206911
Detector/Sample Diluent	101359	212705
Biotinylation Reaction Buffer	101358	214005
Bead Conjugation Buffer	101357	210809
Bead Blocking Buffer	101356	206118
Bead Wash Buffer	101355	206115
Bead Diluent Buffer	100458	210310

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)	1.26 x 10 ⁹
--------------------------	------------------------

Review/Approval

Matt Smart

QC

MAT 31 May 2022

Name

Title

Signature/Date