

Product Number:	101354
Lot Number:	503773
Expiration:	5-Dec-2023

Component	Item Number	Lot Number
488-dyed Singleplex Beads ¹	104006	231817
Detector/Sample Diluent	101359	234016
Biotinylation Reaction Buffer	101358	234206
Bead Conjugation Buffer	101357	233907
Bead Blocking Buffer	101356	234208
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.quoterix.com/HomebrewBeadUpdate>



Bead Count (beads/mL)	1.25 x 10 ⁹
--------------------------	------------------------

Review/Approval

Sonia Khillan *QC Manager* *Sonia Khillan 12/11/22 DEC 2022*

Name

Title

Signature/Date