

| | |
|-----------------|------------|
| Product Number: | 101354 |
| Lot Number: | 503437 |
| Expiration: | 2-Oct-2022 |

| Component | Item Number | Lot Number |
|--|-------------|------------|
| 488-dyed Singleplex Beads ¹ | 104006 | 131609 |
| Detector/Sample Diluent | 101359 | 134925 |
| Biotinylation Reaction Buffer | 101358 | 203014 |
| Bead Conjugation Buffer | 101357 | 134608 |
| Bead Blocking Buffer | 101356 | 127515 |
| Bead Wash Buffer | 101355 | 133218 |
| Bead Diluent Buffer | 100458 | 134609 |

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) | 1.31 x 10 ⁹ |
|--------------------------|------------------------|

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 11 Apr 2022

Name

Title

Signature/Date

Statement: The above product was tested using Quanterix procedures and conforms to Quanterix's specifications.