

| | |
|-----------------|-------------|
| Product Number: | 101354 |
| Lot Number: | 502525 |
| Expiration: | 26-Apr-2021 |

| Component | Item Number | Lot Number |
|-------------------------------|-------------|------------|
| Beads | 103612 | 015313 |
| Detector/Sample Diluent | 101359 | 019711 |
| Biotinylation Reaction Buffer | 101358 | 015108 |
| Bead Conjugation Buffer | 101357 | 011708 |
| Bead Blocking Buffer | 101356 | 019812 |
| Bead Wash Buffer | 101355 | 015106 |
| Bead Diluent Buffer | 100458 | 014512 |

| | |
|--------------------------|----------------------|
| Bead Count (beads/mL) | 3.24x10 ⁹ |
|--------------------------|----------------------|

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 10 Sep 2020
Signature/Date

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