

| | |
|-----------------|-------------|
| Product Number: | 101354 |
| Lot Number: | 502275 |
| Expiration: | 16-Oct-2020 |

| Component | Item Number | Lot Number |
|-------------------------------|-------------|------------|
| Beads | 103207 | 003609 |
| Detector/Sample Diluent | 101359 | 003515 |
| Biotinylation Reaction Buffer | 101358 | 935207 |
| Bead Conjugation Buffer | 101357 | 001506 |
| Bead Blocking Buffer | 101356 | 935209 |
| Bead Wash Buffer | 101355 | 928912 |
| Bead Diluent Buffer | 100458 | 001517 |

| | |
|--------------------------|----------------------|
| Bead Count (beads/mL) | 3.34x10 ⁹ |
|--------------------------|----------------------|

Review/Approval

| | | |
|------------------|--------------------|-------------------------|
| Patrizia Stadler | QC Assoc. Director | <i>P. Stadler</i> |
| _____ Name | _____ Title | _____ Signature/Date |

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