

|                 |            |
|-----------------|------------|
| Product Number: | 102040     |
| Lot Number:     | 502056     |
| Expiration:     | 6-Mar-2020 |

| Component                     | Part Number | Lot Number |
|-------------------------------|-------------|------------|
| Detector/Sample Diluent       | 101359      | 923111     |
| Bead Conjugation Buffer       | 101357      | 909319     |
| Bead Diluent Buffer           | 100458      | 917020     |
| Biotinylation Reaction Buffer | 101358      | 917620     |
| Bead Blocking Buffer          | 101356      | 906502     |
| Bead Wash Buffer              | 101355      | 919220     |

Review/Approval

Patrizia Stadler  
Name

Assoc. Director, QC  
Title

*p. stadler* 09 Oct 2019  
Signature/Date