

| | |
|-----------------|------------|
| Product Number: | 102040 |
| Lot Number: | 503530 |
| Expiration: | 2-Mar-2023 |

| Component | Item Number | Lot Number |
|-------------------------------|-------------|------------|
| Detector/Sample Diluent | 101359 | 214208 |
| Biotinylation Reaction Buffer | 101358 | 209003 |
| Bead Conjugation Buffer | 101357 | 210809 |
| Bead Blocking Buffer | 101356 | 206118 |
| Bead Wash Buffer | 101355 | 206115 |
| Bead Diluent Buffer | 100458 | 210310 |

Review/Approval

Matt Smart
Name

QC

Title

M. Smart 28 Jun 2022
Signature/Date