

| | |
|---------------------|--------------------|
| Product Number: | 104112 |
| Lot Number: | 503883 |
| Paired Reagent Kit: | 503843 |
| Expiration: | 23-Mar-2024 |
| Platform(s): | HD-1 / HD-X & SR-X |

| Component | Part Number | Lot Number | Control Lot Number |
|------------------------|-------------|------------|--------------------|
| Sample Diluent | 104101 | 310104 | 310104A |
| Calibrator Diluent | 104103 | 308217 | N/A |
| Calibrator Concentrate | 104102 | 226601 | N/A |

| | |
|---|-------|
| Calibrator Concentrate Value Assignment (pg/mL) | 47678 |
|---|-------|

Review/Approval

Sonia Khillga QC Sonia Khillga 22 MAY 2023
Name Title Signature/Date