

| | |
|-----------------|-------------|
| Product Number: | 103186 |
| Lot Number: | 503729 |
| Expiration: | 28-Sep-2023 |
| Platform(s): | HD-1 / HD-X |



Data below represents results generated on the Simoa HD-X® Analyzer.

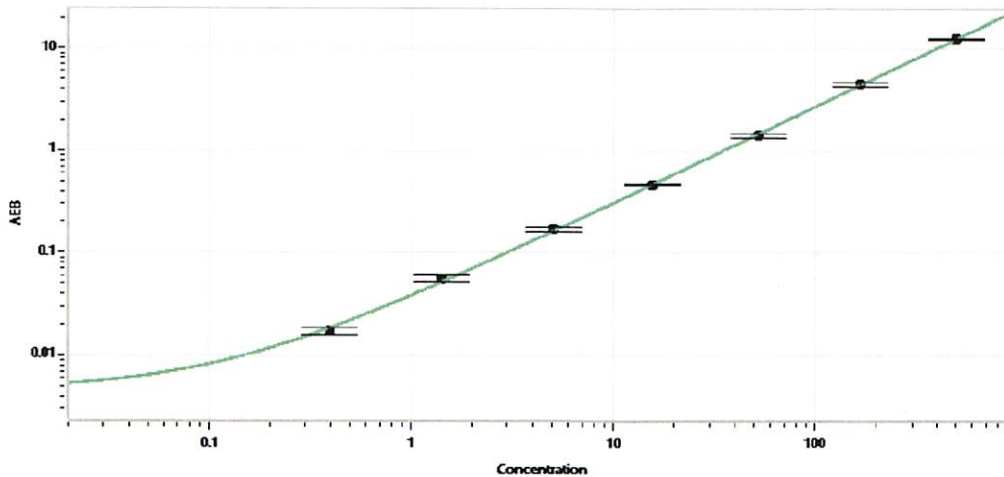
| Component | Part Number | Lot Number |
|------------------|-------------|------------------|
| Bead Reagent | 102246 | 227107 |
| Detector Reagent | 102248 | 227218 |
| SBG Reagent | 102250 | 229301 |
| Sample Diluent | 102252 | 227220 |
| Calibrators | 102255 | 227605 |
| Control 1 | 102256 | 227606 |
| Control 2 | 102257 | 227606 |
| RGP Reagent | 103159 | N/A ¹ |

¹ RGP is not Kit Lot Specific

| Release Materials* | | | |
|--------------------|--------|----------------|---------------------|
| | Lot | Result (pg/mL) | Spec. Range (pg/mL) |
| Control 1 | 227606 | 10.7 | 8.95-13.4 |
| Control 2 | 227606 | 518 | 386-602 |
| Panel 1 | 224402 | 8.69 | 7.80-12.5 |
| Panel 2 | 224402 | 6.80 | 5.71-9.38 |
| Panel 3 | 224402 | 64.8 | 60.3-91.5 |
| Panel 4 | 212511 | 268 | 166-325 |

* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibration Curve



| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| A | 0.000 |
| B | 0.398 |
| C | 1.41 |
| D | 5.04 |
| E | 15.5 |
| F | 52.3 |
| G | 166 |
| H | 493 |

Review/Approval Sonia Khillar QC manager Sonia Khillar / 08 NOV 2022
 Name Title Signature/Date