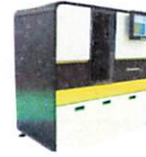


| | |
|------------------------|--------------------|
| Product Number: | 101580 |
| Lot Number: | 504258 |
| Expiration: | 14-Jan-2025 |
| Platform(s): | SR-X and HD-1/HD-X |



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

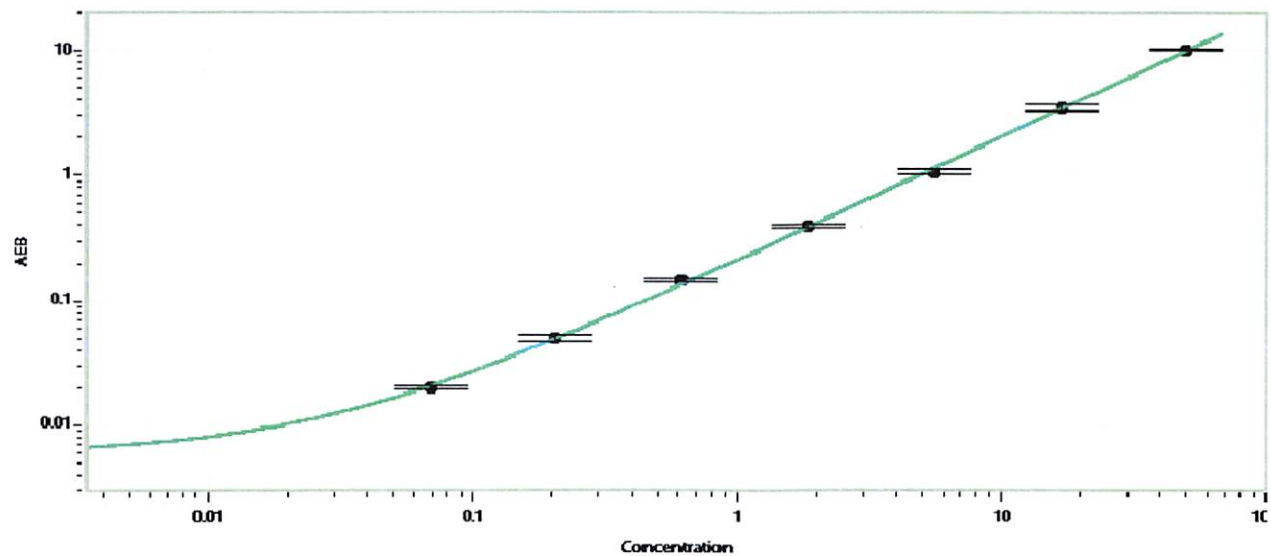
| Component | Part Number | Lot Number | Control Lot Number |
|------------------------|-------------|------------------|--------------------|
| Bead Reagent | 104236 | 319504 | 319504A |
| Detector Reagent | 104327 | 319805 | 319805A |
| SBG Reagent | 104329 | 319806 | 319806A |
| Sample Diluent | 104331 | 319506 | 319506A |
| Calibrator Diluent | 101585 | 319804 | 319804A |
| Calibrator Concentrate | 104332 | 320906 | N/A |
| RGP Reagent | 103159 | N/A ¹ | N/A ¹ |

¹ Reagents are not Kit Lot Specific

| Release Materials* | | | |
|--------------------|--------|----------------|---------------------|
| | Lot | Result (pg/mL) | Spec. Range (pg/mL) |
| Control 1 | 206502 | 3.20 | 2.21 - 3.51 |
| Control 2 | 206502 | 52.4 | 35.6 - 58.8 |
| Panel 2 | 115317 | 2.32 | 1.51 - 3.73 |
| Panel 3 | 116301 | 30.6 | 15.5 - 39.3 |

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

Calibrator Concentrate Value Assignment (pg/mL) 1408



Calibrator Levels (pg/mL)

| | |
|---|-------|
| A | 0.000 |
| B | 0.069 |
| C | 0.206 |
| D | 0.617 |
| E | 1.85 |
| F | 5.56 |
| G | 16.7 |
| H | 50.0 |

Review/Approval

Robert Somogyi
Name

QC Manager
Title

[Signature]
Signature/Date 08/14/2024