

| | |
|------------------------|------------|
| Product Number: | 100988 |
| Lot Number: | 504143 |
| Expiration: | 7-Apr-2025 |
| Platform(s): | HD-1/HD-X |



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

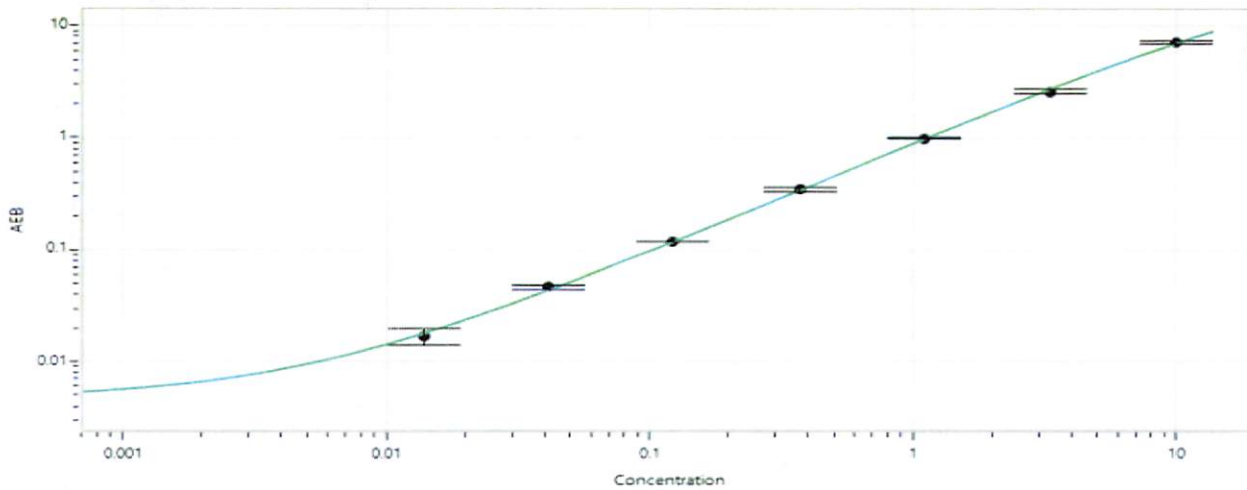
| Component | Part Number | Lot Number |
|------------------------|-------------|------------------|
| Bead Reagent | 100974 | 403110 |
| Detector Reagent | 100976 | 403911 |
| SBG Reagent | 100987 | 403807 |
| Sample Diluent | 103734 | 403813 |
| Calibrator Diluent | 100980 | 407816 |
| Calibrator Concentrate | 100978 | 403912 |
| RGP Reagent | 103159 | N/A ¹ |

¹ Reagents are not Kit Lot Specific

| Release Materials* | | | |
|--------------------|--------|----------------|---------------------|
| | Lot | Result (pg/mL) | Spec. Range (pg/mL) |
| Control 1 | 403912 | 0.093 | 0.075 - 0.125 |
| Control 2 | 403912 | 7.10 | 6.00 - 10.0 |
| Panel 1 | 030001 | 0.195 | 0.105 - 0.234 |
| Panel 2 | 030001 | 0.127 | 0.060 - 0.148 |
| Panel 3 | 030001 | 32.8 | 19.9 - 38.7 |
| Panel 4 | 030001 | 0.202 | 0.111 - 0.215 |

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

| | |
|--|-----|
| Calibrator Concentrate Value Assignment (pg/mL) | 175 |
|--|-----|



| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| A | 0.000 |
| B | 0.014 |
| C | 0.041 |
| D | 0.123 |
| E | 0.370 |
| F | 1.11 |
| G | 3.33 |
| H | 10.0 |

Review/Approval

Cody Carlson
Name

QC Manager
Title

[Signature] 12 APR 2024
Signature/Date