

|                 |                    |
|-----------------|--------------------|
| Product Number: | 104111             |
| Lot Number:     | 503860             |
| Expiration:     | 15-May-2024        |
| Platform(s):    | HD-1 / HD-X & SR-X |



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

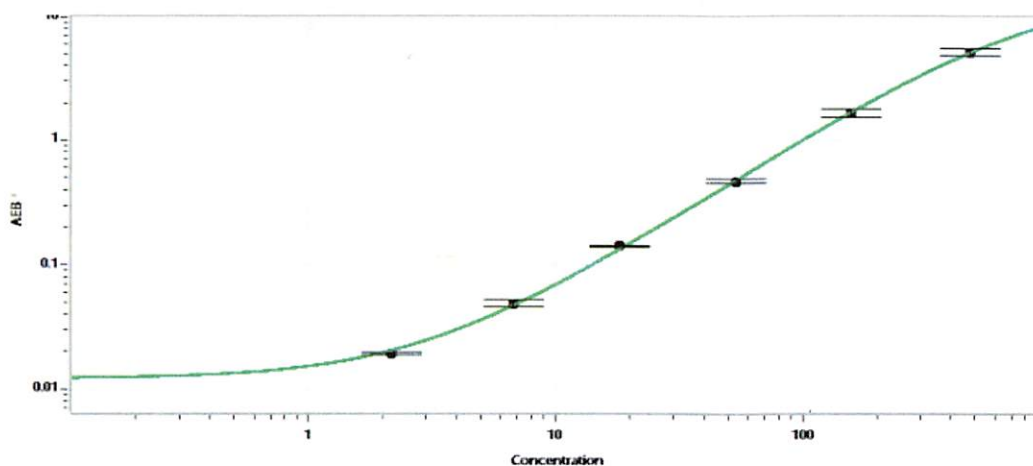
| Component            | Part Number | Lot Number       | Control Lot Number |
|----------------------|-------------|------------------|--------------------|
| Bead Reagent         | 104098      | 313505           | 313505A            |
| Detector Reagent     | 104099      | 313606           | 313606A            |
| SBG Reagent          | 104100      | 313607           | 313607A            |
| Sample Diluent       | 104101      | 313608           | 313608A            |
| Reference Calibrator | 104106      | 315704           | 315704A            |
| Control 1            | 104107      | 315705           | 315705A            |
| Control 2            | 104108      | 315705           | 315705A            |
| RGP Reagent          | 103159      | N/A <sup>1</sup> | N/A <sup>1</sup>   |

<sup>1</sup> RGP is not Kit Lot Specific

| Release Materials* |        |                |                     |
|--------------------|--------|----------------|---------------------|
|                    | Lot    | Result (pg/mL) | Spec. Range (pg/mL) |
| Control 1          | 315705 | 47.6           | 35.2 - 52.8         |
| Control 2          | 315705 | 793            | 697 - 1045          |
| Panel 1            | 301702 | 17.7           | 12.7 - 20.9         |
| Panel 2            | 301702 | 16.9           | 12.3 - 20.9         |
| Panel 3            | 309704 | 458            | 411 - 662           |
| Panel 4            | 301702 | 113            | 57.9 - 125          |

\* 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 | 2.17  |
| C | 6.82  |
| D | 18.2  |
| E | 53.6  |
| F | 156   |
| G | 474   |

Review/Approval

*Sonia Khillan*  
Name

QC

Title

*Sonia Khillan 22 JUN 2023*

Signature/Date