

|                        |             |
|------------------------|-------------|
| Product Number:        | 103769      |
| Lot Number:            | 503181      |
| Expiration:            | 11-Jun-2022 |
| Platform(s):           | HD-X        |
| Product Number (Bulk): | 103845      |
| Lot Number (Bulk):     | N/A         |



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

| Component        | Part Number | Lot Number       |
|------------------|-------------|------------------|
| Bead Reagent     | 103758      | 125701           |
| Detector Reagent | 103760      | 126612           |
| SBG Reagent      | 103762      | 125404           |
| Sample Diluent   | 103763      | 125506           |
| Calibrators      | 103766      | 125803           |
| Control 1        | 103767      | 127903           |
| Control 2        | 103768      | 127903           |
| Control 3        | 103773      | 127903           |
| RGP Reagent      | 103159      | N/A <sup>1</sup> |

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

Release Materials\*

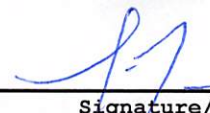
| Sample    | Lot    | Result (ng/mL) | Mean Low limit (ng/mL) | Mean High Limit (ng/mL) |
|-----------|--------|----------------|------------------------|-------------------------|
| Control 1 | 127903 | 380            | 237                    | 493                     |
| Control 2 | 127903 | 3771           | 2918                   | 4377                    |
| Control 3 | 127903 | 18779          | 13634                  | 20451                   |

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

| Sample   | Lot    | Result (ng/mL) | Mean Low limit (ng/mL) | Mean High Limit (ng/mL) |
|----------|--------|----------------|------------------------|-------------------------|
| Panel 1  | 111710 | 0.98           | 0.800                  | 1.29                    |
| Panel 2  | 111710 | 1.94           | 1.51                   | 2.30                    |
| Panel 3  | 111710 | 3.73           | 2.97                   | 4.46                    |
| Panel 4  | 111710 | 6.94           | 5.67                   | 8.50                    |
| Panel 5  | 111710 | 14.4           | 9.74                   | 18.9                    |
| Panel 6  | 111710 | 27.9           | 18.3                   | 36.8                    |
| Panel 8  | 111710 | 6.86           | 4.68                   | 8.48                    |
| Panel 9  | 111710 | 13.6           | 6.86                   | 16.7                    |
| Panel 10 | 111710 | 52.6           | 23.7                   | 67.5                    |
| Panel 11 | 111710 | 0.234          | Reported Result        |                         |
| Panel 12 | 111710 | 0.036          | Reported Result        |                         |
| Panel 13 | 111710 | 0.033          | Reported Result        |                         |
| Panel 14 | 111710 | 0.036          | Reported Result        |                         |
| Panel 15 | 111710 | 0.043          | Reported Result        |                         |
| Panel 16 | 111710 | 0.032          | Reported Result        |                         |
| Panel 17 | 111710 | 0.060          | Reported Result        |                         |
| Panel 18 | 111710 | 0.045          | Reported Result        |                         |
| Panel 19 | 111710 | 0.014          | Reported Result        |                         |
| Panel 20 | 111710 | 0.517          | Reported Result        |                         |

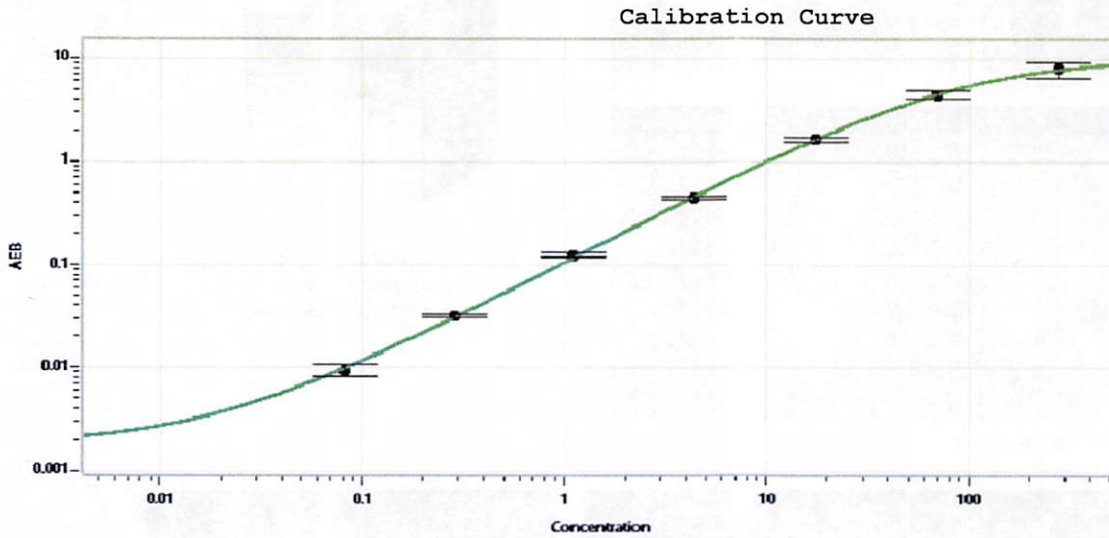
Review/Approval

Wenkai Tsai QC Supervisor  
Name Title

  
Signature/Date

09 Nov 2021





**Calibrator Levels  
(ng/mL)**

|          |       |
|----------|-------|
| <b>A</b> | 0.000 |
| <b>B</b> | 0.082 |
| <b>C</b> | 0.287 |
| <b>D</b> | 1.11  |
| <b>E</b> | 4.39  |
| <b>F</b> | 17.5  |
| <b>G</b> | 69.9  |
| <b>H</b> | 280   |