

| | | |
|-------------------------|--------------------|-------------------|
| Product Number: | 103769 | |
| Lot Number: | 504287 | |
| Expiration Date: | 2-Nov-25 | |
| Platform(s): | HD-X | |
| Component | Part Number | Lot Number |
| Bead Reagent | 103758 | 416301 |
| Detector Reagent | 103760 | 416511 |
| SBG Reagent | 103762 | 419211 |
| Sample Diluent Reagent | 103763 | 416302 |
| Reference Calibrators | 103766 | 416909 |
| Control 1 | 103767 | 416910 |
| Control 2 | 103768 | 416910 |
| Control 3 | 103773 | 416910 |
| RGP Reagent | 103159 | N/A ¹ |



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

¹ Reagents are not Kit Lot Specific

| Release Materials* | | | | |
|--------------------|--------|----------------|------------------------|-------------------------|
| | Lot | Result (ng/mL) | Mean Low Limit (ng/mL) | Mean High Limit (ng/mL) |
| Control 1 | 416910 | 306 | 254 | 459 |
| Control 2 | 416910 | 2671 | 2229 | 4395 |
| Control 3 | 416910 | 12472 | 9364 | 20443 |

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

Review/Approval

Cody Carlson

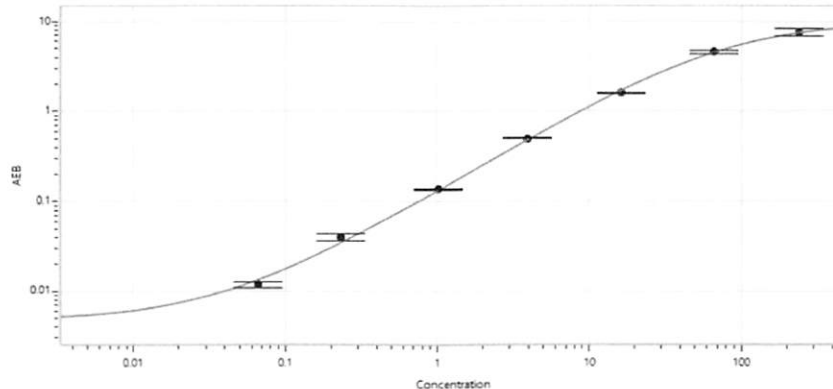
QC Manager

Name

Title

[Handwritten Signature] 09 AUG 2024
Signature/Date

Calibrator Curve



**Calibrator Levels
(pg/mL)**

| | |
|---|-------|
| A | 0.000 |
| B | 0.066 |
| C | 0.233 |
| D | 1.02 |
| E | 3.92 |
| F | 16.3 |
| G | 66.8 |
| H | 244 |