

| | |
|------------------------|-------------|
| Product Number: | 103836 |
| Lot Number: | 504017 |
| Expiration: | 10-Apr-2025 |
| Platform(s): | HD-X |



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

| Component | Part Number | Lot Number |
|---------------------------|-------------|------------------|
| Bead Reagent | 103820 | 328906 |
| Detector Reagent | 103822 | 329003 |
| SBG Reagent | 103824 | 329004 |
| Sample Diluent | 103826 | 328407 |
| Calibrator Diluent | 103829 | 328315 |
| Control Diluent | 104435 | 328316 |
| Calibrator Concentrate | 104443 | 335302 |
| Control 1 Concentrate | 104445 | 335303 |
| Control 2 Concentrate | 104446 | 335304 |
| RGP 2.0 Reagent | 104321 | N/A ¹ |
| Activation Buffer Reagent | 104324 | N/A ¹ |

¹ Reagents are not Kit Lot Specific

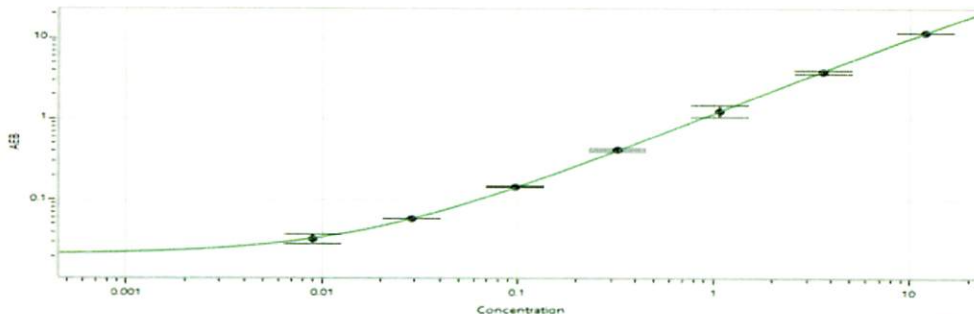
| Release Materials* | | | |
|-----------------------|--------|----------------|---------------------|
| | Lot | Result (pg/mL) | Spec. Range (pg/mL) |
| Control 1 Concentrate | 335303 | 0.192 | 0.074 - 0.251 |
| Control 2 Concentrate | 335304 | 3.80 | 2.15 - 4.00 |
| Panel 1 | 204807 | 0.037 | 0.032 - 0.060 |
| Panel 2 | 204807 | 0.031 | 0.025 - 0.055 |
| Panel 3 | 204807 | 1.42 | 0.932 - 1.73 |
| Panel 4 | 204807 | 2.91 | 1.75 - 3.64 |

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

| | |
|--|------|
| Calibrator Concentrate Reconstitution Volume (uL) | 1045 |
| Control 1 Concentrate Reconstitution Volume (uL) | 1045 |
| Control 2 Concentrate Reconstitution Volume (uL) | 1045 |

| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| A | 0 |
| B | 0.009 |
| C | 0.029 |
| D | 0.098 |
| E | 0.325 |
| F | 1.08 |
| G | 3.60 |
| H | 12.0 |

Calibration Curve



Review/Approval

ROBERT SOMOGYI QC MANAGER

[Handwritten Signature] 03 JAN 2024

Name

Title

Signature/Date