

| | |
|------------------------|-------------|
| Product Number: | 104691 |
| Lot Number: | 504379 |
| Expiration: | 12-Feb-2026 |
| Platform(s): | HD-X |



| Component | Part Number | Lot Number |
|---------------------------|-------------|------------------|
| Bead Reagent | 104752 | 424302 |
| Detector Reagent | 104675 | 424803 |
| SBG Reagent | 104677 | 424804 |
| Plasma Sample Diluent | 104719 | 422524 |
| CSF Sample Diluent | 104720 | 424212 |
| Calibrator Diluent | 104678 | 422624 |
| Control Diluent | 104718 | 422625 |
| Calibrator Concentrate | 104687 | 402944 |
| Control 1 Concentrate | 104689 | 402945 |
| Control 2 Concentrate | 104690 | 402946 |
| RGP 2.0 Reagent | 104321 | N/A ¹ |
| Activation Buffer Reagent | 104324 | N/A ¹ |

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

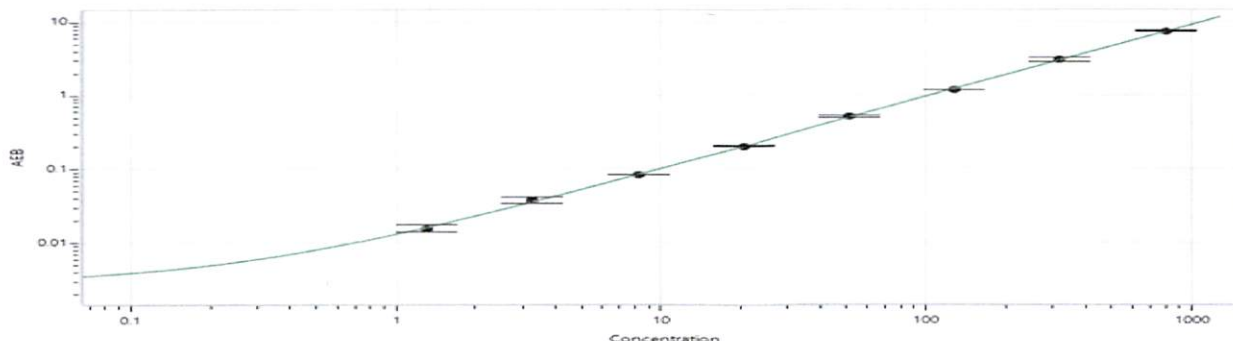
¹ Reagents are not Kit Lot Specific

| Release Materials* | | | |
|-----------------------|--------|----------------|---------------------|
| | Lot | Result (pg/mL) | Spec. Range (pg/mL) |
| Control 1 Concentrate | 402945 | 102 | 77.5 - 144 |
| Control 2 Concentrate | 402946 | 1344 | 978 - 1817 |

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

| | |
|---|------|
| Calibrator Concentrate Reconstitution Volume (uL) | 1037 |
| Control 1 Concentrate Reconstitution Volume (uL) | 1037 |
| Control 2 Concentrate Reconstitution Volume (uL) | 1037 |

Calibration Curve



Calibrator Levels (pg/mL)

| | |
|---|-------|
| A | 0.000 |
| B | 1.31 |
| C | 3.28 |
| D | 8.19 |
| E | 20.5 |
| F | 51.2 |
| G | 128 |
| H | 320 |
| I | 800 |

Review/Approval

Dipesh Patel
Name Title

at 28 Oct 2024
Signature/Date