

|                 |               |
|-----------------|---------------|
| Product Number: | 104364        |
| Lot Number:     | 504674        |
| Expiration:     | 7-May-2027    |
| Platform(s):    | HD-X and SR-X |

| Component                  | Part Number | Lot Number       | Control Lot Number |
|----------------------------|-------------|------------------|--------------------|
| Bead Reagent               | 104352      | 520204           | N/A                |
| Detector Reagent           | 104353      | 506607           | N/A                |
| SBG Reagent                | 104354      | 506610           | 506610A            |
| Sample Diluent Reagent     | 104355      | 506605           | N/A                |
| Calibrator Diluent Reagent | 104358      | 506916           | N/A                |
| Control Diluent Reagent    | 104359      | 506917           | N/A                |
| Calibrator Concentrate     | 104361      | 412713           | N/A                |
| Control 1 Concentrate      | 104362      | 412714           | N/A                |
| Control 2 Concentrate      | 104363      | 412715           | N/A                |
| RGP 2.0 Reagent            | 104321      | N/A <sup>1</sup> | N/A <sup>1</sup>   |
| Activation Buffer Reagent  | 104324      | N/A <sup>1</sup> | N/A <sup>1</sup>   |

<sup>1</sup> Reagents are not Kit Lot Specific



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

| Release Materials* |        |                |                     |
|--------------------|--------|----------------|---------------------|
|                    | Lot    | Result (pg/mL) | Spec. Range (pg/mL) |
| Control 1          | 412714 | 4.41           | 2.98 - 5.53         |
| Control 2          | 412715 | 303            | 242 - 449           |

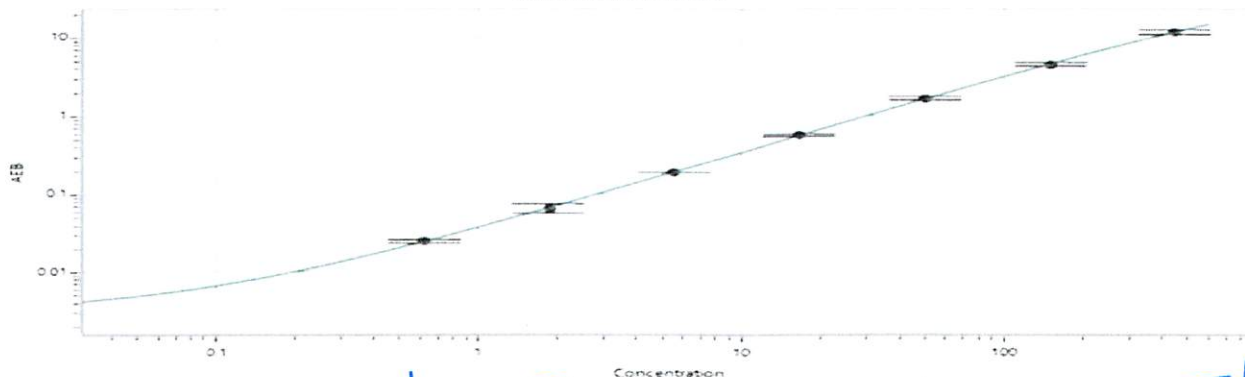
\* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

|                                                   |      |
|---------------------------------------------------|------|
| Calibrator Concentrate Reconstitution Volume (uL) | 1201 |
| Control 1 Concentrate Reconstitution Volume (uL)  | 1201 |
| Control 2 Concentrate Reconstitution Volume (uL)  | 1201 |

**Calibrator Levels (pg/mL)**

|   |       |
|---|-------|
| A | 0.000 |
| B | 0.617 |
| C | 1.85  |
| D | 5.56  |
| E | 16.7  |
| F | 50.0  |
| G | 150   |
| H | 450   |

Calibration Curve



Review/Approval

*Dhruvi Patel*  
Name

*DOC Control II, QA*  
Title

*SR / 13 AUG 2025*  
Signature/Date