

| | |
|-----------------|-------------|
| Product Number: | 105212 |
| Lot Number: | 504425 |
| Expiration: | 15-Apr-2026 |
| Platform(s): | HD-X |



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

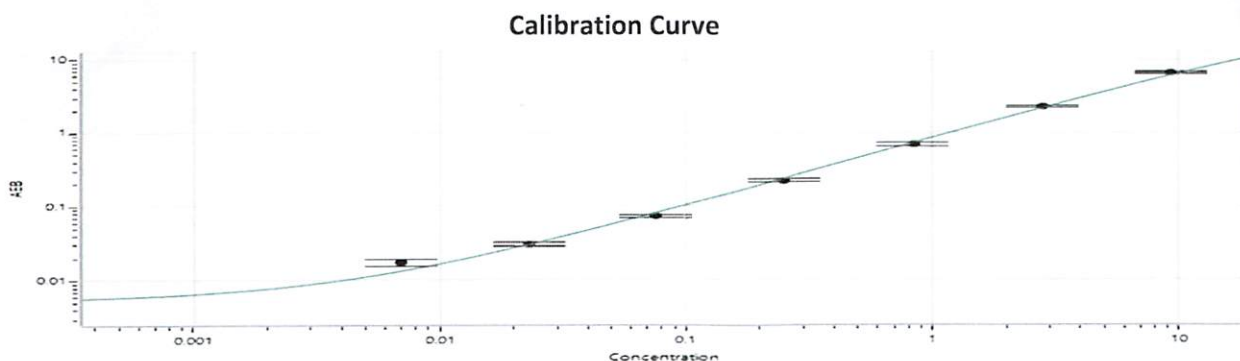
| Component | Part Number | Lot Number |
|---------------------------|-------------|------------------|
| Bead Reagent | 105190 | 429807 |
| Detector Reagent | 105192 | 430202 |
| SBG Reagent | 105194 | 430203 |
| Sample Diluent | 105195 | 429522 |
| Calibrator Diluent | 105197 | 428920 |
| Control Diluent | 105639 | 428921 |
| Calibrator Concentrate | 105208 | 411418 |
| Control 1 Concentrate | 105210 | 411420 |
| Control 2 Concentrate | 105211 | 411421 |
| 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 | 411420 | 0.402 | 0.288 - 0.535 |
| Control 2 Concentrate | 411421 | 23.6 | 16.3 - 30.3 |

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

| | |
|---|------|
| Calibrator Concentrate Reconstitution Volume (uL) | 1000 |
| Control 1 Concentrate Reconstitution Volume (uL) | 1000 |
| Control 2 Concentrate Reconstitution Volume (uL) | 1000 |



| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| A | 0.000 |
| B | 0.007 |
| C | 0.023 |
| D | 0.075 |
| E | 0.252 |
| F | 0.838 |
| G | 2.79 |
| H | 9.29 |

Review/Approval

Dhruvi Patel

DOC Control I, QA

DP 30 Jun 2025

Name

Title

Signature/Date