

|                                |               |
|--------------------------------|---------------|
| Product Number:                | 104671        |
| Lot Number:                    | 504677        |
| Paired Reagent Kit Lot Number: | 504629        |
| Expiration:                    | 22-Jun-2028   |
| Platform(s):                   | HD-X and SR-X |

| Component              | Part Number | Lot Number | Control Lot Number |
|------------------------|-------------|------------|--------------------|
| Calibrator Concentrate | 104666      | 317341     | N/A                |
| Control 1 Concentrate  | 104668      | 317342     | N/A                |
| Control 2 Concentrate  | 104669      | 317343     | 317343A            |



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

|                                                   |      |
|---------------------------------------------------|------|
| Calibrator Concentrate Reconstitution Volume (uL) | 1144 |
| Control 1 Concentrate Reconstitution Volume (uL)  | 1144 |
| Control 2 Concentrate Reconstitution Volume (uL)  | 1144 |

Review/Approval

*Shwami Patel*

Name

*DOC Control II, QA*

Title

*DP/18Jun2025*

Signature/Date