

| | |
|------------------------|------------|
| Product Number: | 100-0217 |
| Lot Number: | 319907 |
| Expiration: | 3-Feb-2024 |
| Platform(s): | SP-X |



Data below represents results generated on the SP-X™ Imaging and Analysis System

| Component | Part Number | Lot Number |
|---------------------------------------|-------------|------------|
| IL-6 Plate | 197-0217 | 319907P |
| Biotinylated Antibody Reagent | 198-0217 | 319907B |
| IL-6 Calibrator | 1861002 | 319285 |
| Sample Diluent 16 | 1863325 | N/A |
| Wash Buffer (12.5X) ¹ | 1863332 | N/A |
| Streptavidin-HRP Reagent ¹ | 1858263 | N/A |
| Super Signal® Luminol ^{1,2} | 1858089 | N/A |
| Super Signal® Peroxide ^{1,2} | 1858090 | N/A |

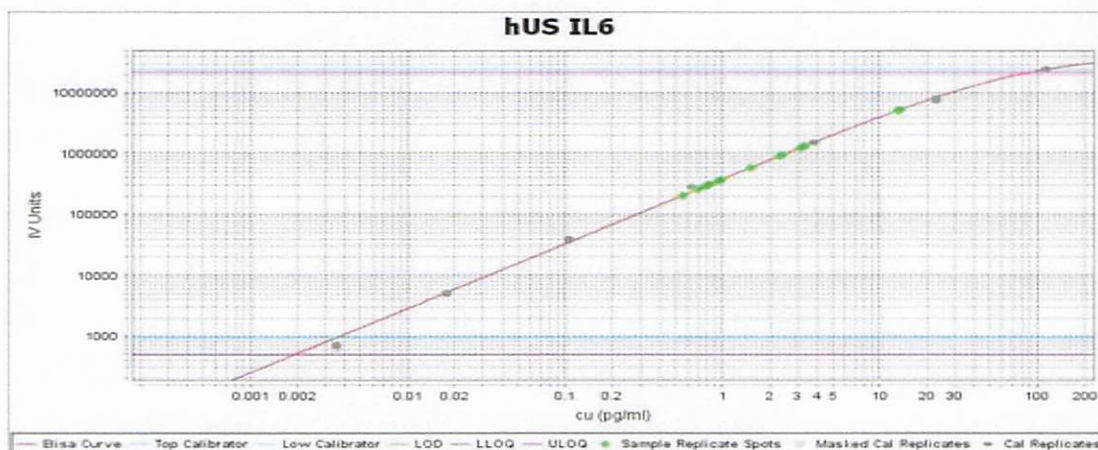
¹ Reagents are not Kit Lot Specific

² SuperSignal® technology is protected by US Patent #6,432,662 and is a registered trademark of ThermoFisher Scientific.

All reagents are specific for use on the SP-X™ platform - refer to the Simoa® IL-6 Developer Kit Instructions (KI-0196).

| Release Materials | | |
|-------------------|----------------|---------------------|
| IL-6 | | |
| | Result (pg/mL) | Spec. Range (pg/mL) |
| Control 2 | 14.2 | 4.53 - 21.5 |
| Control 3 | 3.35 | 1.59 - 5.51 |

IL-6 Calibrator Curve



| Calibrator Levels (pg/mL) | |
|---------------------------|---------|
| 1 | 114 |
| 2 | 22.8 |
| 3 | 3.80 |
| 4 | 0.633 |
| 5 | 0.106 |
| 6 | 0.0176 |
| 7 | 0.00350 |
| 0 | 0.000 |

Review/Approval

Kyle Giroux
Name

Senior Document Control Specialist, QA
Title

Kyle Giroux 26 Feb 2025
Signature/Date