



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 6
Catalog Number: C0072-2
Lot Number: A00U0404
Expiration Date: 31 DEC 2029
Storage Temperature: 2-8°C
Intended Assay Species: Human, NHP

Concentration Assignment:

| Analyte | Amount of Analyte in Calibrator | Unit of Measure | Stock Concentration of Reconstituted Calibrator | Unit of Measure |
|--------------|---------------------------------|-----------------|---|-----------------|
| IL-17A/F | 20625 | pg | 82500 | pg/mL |
| IL-17E/IL-25 | 9062.5 | pg | 36250 | pg/mL |
| IL-17F | 157500 | pg | 630000 | pg/mL |
| IL-21 | 17875 | pg | 71500 | pg/mL |
| IL-22 | 3375 | pg | 13500 | pg/mL |
| IL-23 | 34625 | pg | 138500 | pg/mL |
| IL-27 | 52500 | pg | 210000 | pg/mL |
| IL-29/IFN-λ1 | 10137.5 | pg | 40550 | pg/mL |
| IL-31 | 11800 | pg | 47200 | pg/mL |
| IL-33 | 11387.5 | pg | 45550 | pg/mL |
| TSLP | 9087.5 | pg | 36350 | pg/mL |

Reconstitution Volume: 250 µL

To calculate the **Calibrator Standard 1 (top of curve) concentration**, please refer to the dilution-factor used in the **Prepare Calibrator Standards** section of the respective product insert. Divide the stock concentration of the reconstituted calibrator by this dilution factor to generate the working calibrator standard 1 concentration.

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

| | Name | Function | Signature | Date |
|------------------------|--------------|----------|-----------|-------------|
| Review/Approval | Karen Hamila | Quality | | 04 MAR 2025 |