



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 9
Catalog Number: C0090-2
Lot Number: A00U0351
Expiration Date: 30 NOV 2028
Storage Temperature: 2-8°C
Intended Assay Species: Human

Concentration Assignment:

| Analyte | Amount of Analyte in Calibrator | Unit of Measure | Stock Concentration of Reconstituted Calibrator | Unit of Measure |
|----------------|---------------------------------|-----------------|---|-----------------|
| EPO | 26250 | pg | 105000 | pg/mL |
| FLT3L | 7750 | pg | 31000 | pg/mL |
| IFN- β | 122500 | pg | 490000 | pg/mL |
| IL-3 | 19000 | pg | 76000 | pg/mL |
| IL-9 | 1662.5 | pg | 6650 | pg/mL |
| IL-17B | 3425 | pg | 13700 | pg/mL |
| IL-17C | 24250 | pg | 97000 | pg/mL |
| IL-17D | 31500 | pg | 126000 | pg/mL |
| IL-1RA | 6225 | pg | 24900 | pg/mL |
| IL-2R α | 71000 | pg | 284000 | pg/mL |

Reconstitution Volume: 250 μ L

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 Hamilla | Quality | | 12 JAN 2024 |