



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 9

Catalog Number: C0090-2

Lot Number: A00U0258

Expiration Date: 31 MAY 2027

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 | 25625 | pg | 102500 | pg/mL |
| FLT3L | 7238 | pg | 28950 | pg/mL |
| IFN- β | 135000 | pg | 540000 | pg/mL |
| IL-3 | 18875 | pg | 75500 | pg/mL |
| IL-9 | 1638 | pg | 6550 | pg/mL |
| IL-17B | 4100 | pg | 16400 | pg/mL |
| IL-17C | 19375 | pg | 77500 | pg/mL |
| IL-17D | 33250 | pg | 133000 | pg/mL |
| IL-IRA | 5875 | pg | 23500 | pg/mL |
| IL-2R α | 69625 | pg | 278500 | 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 | | 15 JUL 2022 |