



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 9

**Catalog Number:** C0090-2

**Lot Number:** A00U0230

**Expiration Date:** 31 JAN 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	24000	pg	96000	pg/mL
FLT3L	6913	pg	27650	pg/mL
IFN- $\beta$	130000	pg	520000	pg/mL
IL-3	18875	pg	75500	pg/mL
IL-9	1488	pg	5950	pg/mL
IL-17B	4663	pg	18650	pg/mL
IL-17C	19750	pg	79000	pg/mL
IL-17D	38625	pg	154500	pg/mL
IL-IRA	5738	pg	22950	pg/mL
IL-2R $\alpha$	58250	pg	233000	pg/mL

**Reconstitution Volume:** 250  $\mu$ 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
<b>Review/Approval</b>	Karen Hamilla	Quality		23 MAR 2022