



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 9  
**Catalog Number:** C0090-2  
**Lot Number:** A00U0371  
**Expiration Date:** 31 MAR 2029  
**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	26750	pg	107000	pg/mL
FLT3L	7862.5	pg	31450	pg/mL
IFN- $\beta$	12625	pg	50500	pg/mL
IL-3	19750	pg	79000	pg/mL
IL-9	1712.5	pg	6850	pg/mL
IL-17B	3337.5	pg	13350	pg/mL
IL-17C	24875	pg	99500	pg/mL
IL-17D	37750	pg	151000	pg/mL
IL-1RA	5687.5	pg	22750	pg/mL
IL-2R $\alpha$	72375	pg	289500	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		15 MAY 2024