



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 9  
**Catalog Number:** C0090-2  
**Lot Number:** A00U0210  
**Expiration Date:** 30 APR 2026  
**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	24250	pg	97000	pg/mL
FLT3L	7500	pg	30000	pg/mL
IFN-β	126250	pg	505000	pg/mL
IL-3	20625	pg	82500	pg/mL
IL-9	1700	pg	6800	pg/mL
IL-17B	4463	pg	17850	pg/mL
IL-17C	28750	pg	115000	pg/mL
IL-17D	71250	pg	285000	pg/mL
IL-1RA	6488	pg	25950	pg/mL
IL-2Rα	68125	pg	272500	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
<b>Review/Approval</b>	Karen Hamilla	Quality		01 JUL 2021