



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 9

**Catalog Number:** C0090-2

**Lot Number:** A00U0193

**Expiration Date:** 31 JAN 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	24875	pg	99500	pg/mL
FLT3L	7813	pg	31250	pg/mL
IFN- $\beta$	136250	pg	545000	pg/mL
IL-3	21500	pg	86000	pg/mL
IL-9	1713	pg	6850	pg/mL
IL-17B	3450	pg	13800	pg/mL
IL-17C	27125	pg	108500	pg/mL
IL-17D	44625	pg	178500	pg/mL
IL-1RA	5825	pg	23300	pg/mL
IL-2R $\alpha$	67000	pg	268000	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		31 MAR 2021