



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 9

**Catalog Number:** C0090-2

**Lot Number:** A00U0155

**Expiration Date:** 30 APR 2025

**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	24125	pg	96500	pg/mL
FLT3L	7525	pg	30100	pg/mL
IFN- $\beta$	135000	pg	540000	pg/mL
IL-3	21250	pg	85000	pg/mL
IL-9	1563	pg	6250	pg/mL
IL-17B	4975	pg	19900	pg/mL
IL-17C	25375	pg	101500	pg/mL
IL-17D	62250	pg	249000	pg/mL
IL-1RA	6650	pg	26600	pg/mL
IL-2R $\alpha$	66375	pg	265500	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		27 APR 2020