



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 9

Catalog Number: C0090-2

Lot Number: A00U0221

Expiration Date: 30 JUN 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	24000	pg	96000	pg/mL
FLT3L	7588	pg	30350	pg/mL
IFN- β	128750	pg	515000	pg/mL
IL-3	20000	pg	80000	pg/mL
IL-9	1650	pg	6600	pg/mL
IL-17B	4625	pg	18500	pg/mL
IL-17C	33000	pg	132000	pg/mL
IL-17D	85000	pg	340000	pg/mL
IL-1RA	5925	pg	23700	pg/mL
IL-2R α	68750	pg	275000	pg/mL

Reconstitution Volume: 250 μ 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
Review/Approval	Xiomara Talavera	Quality		03 SEP 2021