



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 9

Catalog Number: C0090-2

Lot Number: A00U0165

Expiration Date: 31 AUG 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	25000	pg	100000	pg/mL
FLT3L	7288	pg	29150	pg/mL
IFN- β	131250	pg	525000	pg/mL
IL-3	20750	pg	83000	pg/mL
IL-9	1688	pg	6750	pg/mL
IL-17B	3650	pg	14600	pg/mL
IL-17C	24375	pg	97500	pg/mL
IL-17D	48375	pg	193500	pg/mL
IL-1RA	6000	pg	24000	pg/mL
IL-2R α	66000	pg	264000	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	Karen Hamilla	Quality		04 SEP 2020