



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 9
Catalog Number: C0090-2
Lot Number: A00U0413
Expiration Date: 30 SEP 2030
Storage Temperature: 2-8°C
Intended Assay Species: Human, NHP


Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
EPO	20625	pg	82500	pg/mL
FLT3L	6212.5	pg	24850	pg/mL
IFN- β	12625	pg	50500	pg/mL
IL-3	18250	pg	73000	pg/mL
IL-9	1537.5	pg	6150	pg/mL
IL-17B	4325	pg	17300	pg/mL
IL-17C	20875	pg	83500	pg/mL
IL-17D	45000	pg	180000	pg/mL
IL-1RA	5887.5	pg	23550	pg/mL
IL-2R α	64750	pg	259000	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 DEC 2025