



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 9
Catalog Number: C0090-2
Lot Number: A00U0469
Expiration Date: 31 MAR 2031
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	21250	pg	85000	pg/mL
FLT3L	6325	pg	25300	pg/mL
IFN- β	13375	pg	53500	pg/mL
IL-3	18775	pg	75100	pg/mL
IL-9	1582.5	pg	6330	pg/mL
IL-17B	4825	pg	19300	pg/mL
IL-17C	22025	pg	88100	pg/mL
IL-17D	46000	pg	184000	pg/mL
IL-1RA	5875	pg	23500	pg/mL
IL-2R α	66250	pg	265000	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		30 APR 2026