



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 9
Catalog Number: C0090-2
Lot Number: A00U0405
Expiration Date: 31 JAN 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	21125	pg	84500	pg/mL
FLT3L	7700	pg	30800	pg/mL
IFN- β	12875	pg	51500	pg/mL
IL-3	18250	pg	73000	pg/mL
IL-9	1562.5	pg	6250	pg/mL
IL-17B	2850	pg	11400	pg/mL
IL-17C	20875	pg	83500	pg/mL
IL-17D	39250	pg	157000	pg/mL
IL-1RA	5712.5	pg	22850	pg/mL
IL-2R α	66125	pg	264500	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		10 MAR 2025