



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 9

Catalog Number: C0090-2

Lot Number: A00U0277

Expiration Date: 30 JUN 2027

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	25375	pg	101500	pg/mL
FLT3L	7363	pg	29450	pg/mL
IFN- β	132500	pg	530000	pg/mL
IL-3	18625	pg	74500	pg/mL
IL-9	1663	pg	6650	pg/mL
IL-17B	4275	pg	17100	pg/mL
IL-17C	20250	pg	81000	pg/mL
IL-17D	27500	pg	110000	pg/mL
IL-1RA	5900	pg	23600	pg/mL
IL-2R α	70750	pg	283000	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		14 SEP 2022