



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 9

Catalog Number: C0090-2

Lot Number: A00U0113

Expiration Date: 31 MAY 2024

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	25 500	pg	102 000	pg/mL
FLT3L	7 700	pg	30 800	pg/mL
IFN- β	122 125	pg	488 500	pg/mL
IL-3	17 500	pg	70 000	pg/mL
IL-9	1 675	pg	6 700	pg/mL
IL-17B	3 438	pg	13 750	pg/mL
IL-17C	30 000	pg	120 000	pg/mL
IL-17D	49 000	pg	196 000	pg/mL
IL-1RA	5 813	pg	23 250	pg/mL
IL-2R α	66 625	pg	266 500	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 2019