



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 9

Catalog Number: C0090-2

Lot Number: A00U0144

Expiration Date: 28 FEB 2025

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	26750	pg	107000	pg/mL
FLT3L	7013	pg	28050	pg/mL
IFN- β	122875	pg	491500	pg/mL
IL-3	19500	pg	78000	pg/mL
IL-9	1625	pg	6500	pg/mL
IL-17B	3213	pg	12850	pg/mL
IL-17C	23500	pg	94000	pg/mL
IL-17D	49500	pg	198000	pg/mL
IL-1RA	6863	pg	27450	pg/mL
IL-2R α	69000	pg	276000	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		17 MAR 2020