



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 9
Catalog Number: C0090-2
Lot Number: A00U0302
Expiration Date: 31 DEC 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	26875	pg	107500	pg/mL
FLT3L	7875	pg	31500	pg/mL
IFN- β	135000	pg	540000	pg/mL
IL-3	23375	pg	93500	pg/mL
IL-9	1700	pg	6800	pg/mL
IL-17B	4450	pg	17800	pg/mL
IL-17C	25625	pg	102500	pg/mL
IL-17D	36500	pg	146000	pg/mL
IL-1RA	6012.5	pg	24050	pg/mL
IL-2R α	74250	pg	297000	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 was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		26 JAN 2023