



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 9

Catalog Number: C0090-2

Lot Number: A00U0117

Expiration Date: 31 OCT 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	27250	pg	109000	pg/mL
FLT3L	7900	pg	31600	pg/mL
IFN- β	128750	pg	515000	pg/mL
IL-3	20250	pg	81000	pg/mL
IL-9	1663	pg	6650	pg/mL
IL-17B	4488	pg	17950	pg/mL
IL-17C	27125	pg	108500	pg/mL
IL-17D	43250	pg	173000	pg/mL
IL-1RA	5888	pg	23550	pg/mL
IL-2R α	71125	pg	284500	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		08 NOV 2019