



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 9  
**Catalog Number:** C0090-2  
**Lot Number:** A00U0127  
**Expiration Date:** 30 SEP 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	24 750	pg	99 000	pg/mL
FLT3L	8 225	pg	32 900	pg/mL
IFN- $\beta$	130 000	pg	520 000	pg/mL
IL-3	18 750	pg	75 000	pg/mL
IL-9	1 600	pg	6 400	pg/mL
IL-17B	5 563	pg	22 250	pg/mL
IL-17C	27 375	pg	109 500	pg/mL
IL-17D	59 375	pg	237 500	pg/mL
IL-1RA	6 025	pg	24 100	pg/mL
IL-2R $\alpha$	63 375	pg	253 500	pg/mL

**Reconstitution Volume:** 250  $\mu$ 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
<b>Review/Approval</b>	Xiomara Talavera	Quality		27 SEP 2019