



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 9

**Catalog Number:** C0090-2

**Lot Number:** A00U0103

**Expiration Date:** 30 APR 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	25 625	pg	102 500	pg/mL
FLT3L	8 025	pg	32 100	pg/mL
IFN- $\beta$	114 375	pg	457 500	pg/mL
IL-3	18 875	pg	75 500	pg/mL
IL-9	1 613	pg	6 450	pg/mL
IL-17B	4 313	pg	17 250	pg/mL
IL-17C	27 750	pg	111 000	pg/mL
IL-17D	57 500	pg	230 000	pg/mL
IL-1RA	6 050	pg	24 200	pg/mL
IL-2R $\alpha$	65 000	pg	260 000	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>	Dawn Stollar	Quality		11 APR 2019