



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 1  
**Catalog Number:** C0060-2  
**Lot Number:** A00U0140  
**Expiration Date:** 31 JAN 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
IFN- $\gamma$	41625	pg	166500	pg/mL
IL-1 $\beta$	5625	pg	22500	pg/mL
IL-2	2688	pg	10750	pg/mL
IL-4	3250	pg	13000	pg/mL
IL-6	2875	pg	11500	pg/mL
IL-8	2750	pg	11000	pg/mL
IL-10	4925	pg	19700	pg/mL
IL-12p70	8625	pg	34500	pg/mL
IL-13	3350	pg	13400	pg/mL
TNF- $\alpha$	5600	pg	22400	pg/mL
IL-5	5488	pg	21950	pg/mL
GM-CSF	15250	pg	61000	pg/mL
IL-17A	28500	pg	114000	pg/mL
VEGF-A	9650	pg	38600	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>	Karen Hamilla	Quality		26 FEB 2020