



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 1  
**Catalog Number:** C0060-2  
**Lot Number:** A00U0109  
**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
IFN- $\gamma$	31 125	pg	124 500	pg/mL
IL-1 $\beta$	4 400	pg	17 600	pg/mL
IL-2	2 525	pg	10 100	pg/mL
IL-4	3 050	pg	12 200	pg/mL
IL-6	2 400	pg	9 600	pg/mL
IL-8	2 475	pg	9 900	pg/mL
IL-10	4 438	pg	17 750	pg/mL
IL-12p70	6 850	pg	27 400	pg/mL
IL-13	2 775	pg	11 100	pg/mL
TNF- $\alpha$	4 575	pg	18 300	pg/mL
IL-5	4 825	pg	19 300	pg/mL
GM-CSF	13 375	pg	53 500	pg/mL
IL-17A	24 875	pg	99 500	pg/mL
VEGF-A	8 263	pg	33 050	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		02 OCT 2019