



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 3
Catalog Number: C0062-2
Lot Number: A00U0342
Expiration Date: 31 AUG 2028
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
IL-1 α	6912.5	pg	27650	pg/mL
IL-7	6200	pg	24800	pg/mL
IL-12/IL-23p40	32375	pg	129500	pg/mL
IL-15	3725	pg	14900	pg/mL
IL-16	27500	pg	110000	pg/mL
TNF- β	4150	pg	16600	pg/mL
G-CSF	11975	pg	47900	pg/mL
IFN- α 2a	49125	pg	196500	pg/mL
IL-18	16750	pg	67000	pg/mL
TPO	47875	pg	191500	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 was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		12 SEP 2023