



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 3
Catalog Number: C0062-2
Lot Number: A00U0305
Expiration Date: 31 MAR 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	6975	pg	27900	pg/mL
IL-7	8700	pg	34800	pg/mL
IL-12/IL-23p40	32875	pg	131500	pg/mL
IL-15	3750	pg	15000	pg/mL
IL-16	26250	pg	105000	pg/mL
TNF- β	3837.5	pg	15350	pg/mL
G-CSF	13375	pg	53500	pg/mL
IFN- α 2a	45000	pg	180000	pg/mL
IL-18	14625	pg	58500	pg/mL
TPO	51250	pg	205000	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		09 MAY 2023