



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 3  
**Catalog Number:** C0062-2  
**Lot Number:** A00U0338  
**Expiration Date:** 31 JUL 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 $\alpha$	7475	pg	29900	pg/mL
IL-7	8350	pg	33400	pg/mL
IL-12/IL-23p40	32750	pg	131000	pg/mL
IL-15	3550	pg	14200	pg/mL
IL-16	25625	pg	102500	pg/mL
TNF- $\beta$	3725	pg	14900	pg/mL
G-CSF	12750	pg	51000	pg/mL
IFN- $\alpha$ 2a	44750	pg	179000	pg/mL
IL-18	16125	pg	64500	pg/mL
TPO	46250	pg	185000	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 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
<b>Review/Approval</b>	Karen Hamilla	Quality		07 SEP 2023