



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 5
Catalog Number: C0065-2
Lot Number: A00U0326
Expiration Date: 30 APR 2028
Storage Temperature: 2-8°C
Intended Assay Species: Mouse

Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	3587.5	pg	14350	pg/mL
IL-1 β	16875	pg	67500	pg/mL
IL-2	11825	pg	47300	pg/mL
IL-4	12875	pg	51500	pg/mL
IL-5	3750	pg	15000	pg/mL
IL-6	20375	pg	81500	pg/mL
KC/GRO	2612.5	pg	10450	pg/mL
IL-10	28500	pg	114000	pg/mL
IL-12p70	137500	pg	550000	pg/mL
TNF- α	8700	pg	34800	pg/mL
GM-CSF	1362.5	pg	5450	pg/mL
IL-13	23000	pg	92000	pg/mL
EPO	17500	pg	70000	pg/mL
VEGF-A	13125	pg	52500	pg/mL

Reconstitution Volume: 250 μ L 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	Xiomara Talavera	Quality		25 MAY 2023