



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 5
Catalog Number: C0065-2
Lot Number: A00U0359
Expiration Date: 31 MAR 2029
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	3650	pg	14600	pg/mL
IL-1 β	18375	pg	73500	pg/mL
IL-2	11825	pg	47300	pg/mL
IL-4	11462.5	pg	45850	pg/mL
IL-5	3687.5	pg	14750	pg/mL
IL-6	21500	pg	86000	pg/mL
KC/GRO	2600	pg	10400	pg/mL
IL-10	27625	pg	110500	pg/mL
IL-12p70	127500	pg	510000	pg/mL
TNF- α	7675	pg	30700	pg/mL
GM-CSF	1362.5	pg	5450	pg/mL
IL-13	27375	pg	109500	pg/mL
EPO	16500	pg	66000	pg/mL
VEGF-A	11625	pg	46500	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 is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		09 MAY 2024