



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 5
Catalog Number: C0065-2
Lot Number: A00U0452
Expiration Date: 30 NOV 2030
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	3900	pg	15600	pg/mL
IL-1 β	18425	pg	73700	pg/mL
IL-2	11925	pg	47700	pg/mL
IL-4	12325	pg	49300	pg/mL
IL-5	3625	pg	14500	pg/mL
IL-6	20475	pg	81900	pg/mL
KC/GRO	2312.5	pg	9250	pg/mL
IL-10	30500	pg	122000	pg/mL
IL-12p70	133500	pg	534000	pg/mL
TNF- α	8425	pg	33700	pg/mL
GM-CSF	1232.5	pg	4930	pg/mL
IL-13	23625	pg	94500	pg/mL
IL-13	23625	pg	94500	pg/mL
VEGF-A	12300	pg	49200	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		03 FEB 2026