



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 5

Catalog Number: C0065-2

Lot Number: A00U0119

Expiration Date: 31 JUL 2024

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	3 675	pg	14 700	pg/mL
IL-1 β	12 875	pg	51 500	pg/mL
IL-2	12 750	pg	51 000	pg/mL
IL-4	13 000	pg	52 000	pg/mL
IL-5	3 450	pg	13 800	pg/mL
IL-6	22 375	pg	89 500	pg/mL
KC	2 650	pg	10 600	pg/mL
IL-10	28 500	pg	114 000	pg/mL
IL-12p70	118 125	pg	472 500	pg/mL
TNF- α	8 463	pg	33 850	pg/mL
GM-CSF	1 313	pg	5 250	pg/mL
IL-13	22 750	pg	91 000	pg/mL
EPO	15 750	pg	63 000	pg/mL
VEGF-A	13 250	pg	53 000	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		22 AUG 2019