



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 2
Catalog Number: C0061-2
Lot Number: A00U0369
Expiration Date: 30 APR 2029
Storage Temperature: 2-8°C
Intended Assay Species:

Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Eotaxin	6262.5	pg	25050	pg/mL
Eotaxin-3	23750	pg	95000	pg/mL
IP-10	13500	pg	54000	pg/mL
MCP-1	6750	pg	27000	pg/mL
MCP-4	4812.5	pg	19250	pg/mL
MDC	24000	pg	96000	pg/mL
MIP-1 α	8937.5	pg	35750	pg/mL
MIP-1 β	2187.5	pg	8750	pg/mL
TARC/CCL17	3162.5	pg	12650	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	Sophie Brainerd	Quality		10 JUN 2024