



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 2
Catalog Number: C0061-2
Lot Number: A00U0301
Expiration Date: 30 NOV 2027
Storage Temperature: 2-8°C
Intended Assay Species: Human

Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Eotaxin	9837.5	pg	39350	pg/mL
Eotaxin-3	28875	pg	115500	pg/mL
IP-10	14125	pg	56500	pg/mL
MCP-1	7912.5	pg	31650	pg/mL
MCP-4	4750	pg	19000	pg/mL
MDC	28125	pg	112500	pg/mL
MIP-1 α	7187.5	pg	28750	pg/mL
MIP-1 β	2487.5	pg	9950	pg/mL
TARC/CCL17	3600	pg	14400	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 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	Karen Hamilla	Quality		27 DEC 2022



Page 2 of 2

CERTIFICATE OF ANALYSIS

COA-17609-A

QA-FM-184-F