



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 2
Catalog Number: C0061-2
Lot Number: A00U0337
Expiration Date: 31 JUL 2028
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	7287.5	pg	29150	pg/mL
Eotaxin-3	23250	pg	93000	pg/mL
IP-10	11400	pg	45600	pg/mL
MCP-1	7375	pg	29500	pg/mL
MCP-4	4537.5	pg	18150	pg/mL
MDC	25500	pg	102000	pg/mL
MIP-1 α	7162.5	pg	28650	pg/mL
MIP-1 β	2850	pg	11400	pg/mL
TARC/CCL17	3562.5	pg	14250	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		05 SEP 2023