



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 2

Catalog Number: C0061-2

Lot Number: A00U0275

Expiration Date: 31 JUL 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	11813	pg	47250	pg/mL
Eotaxin-3	29750	pg	119000	pg/mL
IP-10	15750	pg	63000	pg/mL
MCP-1	9538	pg	38150	pg/mL
MCP-4	6138	pg	24550	pg/mL
MDC	32500	pg	130000	pg/mL
MIP-1 α	9263	pg	37050	pg/mL
MIP-1 β	3250	pg	13000	pg/mL
TARC	4600	pg	18400	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		02 SEP 2022