



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 2

Catalog Number: C0061-2

Lot Number: A00U0246

Expiration Date: 31 DEC 2026

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	8863	pg	35450	pg/mL
Eotaxin-3	23125	pg	92500	pg/mL
IP-10	11513	pg	46050	pg/mL
MCP-1	6813	pg	27250	pg/mL
MCP-4	4463	pg	17850	pg/mL
MDC	22125	pg	88500	pg/mL
MIP-1 α	6975	pg	27900	pg/mL
MIP-1 β	2238	pg	8950	pg/mL
TARC	3350	pg	13400	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		11 APR 2022