



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 2

Catalog Number: C0061-2

Lot Number: A00U0256

Expiration Date: 31 MAY 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	9775	pg	39100	pg/mL
Eotaxin-3	26250	pg	105000	pg/mL
IP-10	12875	pg	51500	pg/mL
MCP-1	8000	pg	32000	pg/mL
MCP-4	4963	pg	19850	pg/mL
MDC	25625	pg	102500	pg/mL
MIP-1 α	7188	pg	28750	pg/mL
MIP-1 β	2538	pg	10150	pg/mL
TARC	3838	pg	15350	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	Xiomara Talavera	Quality		01 JUL 2022