



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 2

Catalog Number: C0061 -2

Lot Number: A00U0287

Expiration Date: 30 SEP 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	8313	pg	33250	pg/mL
Eotaxin-3	28500	pg	114000	pg/mL
IP-10	13875	pg	55500	pg/mL
MCP-1	7625	pg	30500	pg/mL
MCP-4	4775	pg	19100	pg/mL
MDC	29125	pg	116500	pg/mL
MIP-1 α	7288	pg	29150	pg/mL
MIP-1 β	3013	pg	12050	pg/mL
TARC	3850	pg	15400	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 NOV 2022