



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 2

Catalog Number: C0061-2

Lot Number: A00U0202

Expiration Date: 28 FEB 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	7875	pg	31500	pg/mL
Eotaxin-3	24250	pg	97000	pg/mL
IP-10	11925	pg	47700	pg/mL
MCP-1	6813	pg	27250	pg/mL
MCP-4	4850	pg	19400	pg/mL
MDC	22875	pg	91500	pg/mL
MIP-1 α	7125	pg	28500	pg/mL
MIP-1 β	2275	pg	9100	pg/mL
TARC	3563	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 is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		15 APR 2021