



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 2

Catalog Number: C0061-2

Lot Number: A00U0142

Expiration Date: 31 JAN 2025

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	7825	pg	31300	pg/mL
Eotaxin-3	24375	pg	97500	pg/mL
IP-10	14625	pg	58500	pg/mL
MCP-1	6800	pg	27200	pg/mL
MCP-4	5225	pg	20900	pg/mL
MDC	29500	pg	118000	pg/mL
MIP-1 α	8988	pg	35950	pg/mL
MIP-1 β	2375	pg	9500	pg/mL
TARC	3488	pg	13950	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		05 MAR 2020