



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 2
Catalog Number: C0061-2
Lot Number: A00U0153
Expiration Date: 30 JUN 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	7188	pg	28750	pg/mL
Eotaxin-3	23750	pg	95000	pg/mL
IP-10	13625	pg	54500	pg/mL
MCP-1	6738	pg	26950	pg/mL
MCP-4	4800	pg	19200	pg/mL
MDC	21500	pg	86000	pg/mL
MIP-1 α	8400	pg	33600	pg/mL
MIP-1 β	2213	pg	8850	pg/mL
TARC	3663	pg	14650	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		29 JUL 2020