



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 2
Catalog Number: C0061-2
Lot Number: A00U0123
Expiration Date: 31 AUG 2024
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	7 413	pg	29 650	pg/mL
Eotaxin-3	24 875	pg	99 500	pg/mL
IP-10	12 750	pg	51 000	pg/mL
MCP-1	6 750	pg	27 000	pg/mL
MCP-4	5 238	pg	20 950	pg/mL
MDC	28 000	pg	112 000	pg/mL
MIP-1 α	8 375	pg	33 500	pg/mL
MIP-1 β	2 288	pg	9 150	pg/mL
TARC	3 563	pg	14 250	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		18 SEP 2019