



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 2

Catalog Number: C0061-2

Lot Number: A00U0130

Expiration Date: 30 NOV 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	7238	pg	28950	pg/mL
Eotaxin-3	24125	pg	96500	pg/mL
IP-10	12875	pg	51500	pg/mL
MCP-1	7113	pg	28450	pg/mL
MCP-4	5125	pg	20500	pg/mL
MDC	27125	pg	108500	pg/mL
MIP-1 α	8125	pg	32500	pg/mL
MIP-1 β	2275	pg	9100	pg/mL
TARC	3688	pg	14750	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		09 JAN 2020