



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 12

Catalog Number: C0092-2

Lot Number: A00U0212

Expiration Date: 30 APR 2026

Storage Temperature: 2-8°C

Intended Assay Species: Mouse

Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IL-9	12488	pg	49950	pg/mL
IL-17A/F	12875	pg	51500	pg/mL
IP-10	7525	pg	30100	pg/mL
MCP-1	1488	pg	5950	pg/mL
MIP-1 α	2200	pg	8800	pg/mL
MIP-1 β	36125	pg	144500	pg/mL
MIP-2	3063	pg	12250	pg/mL
MIP-3 α	3188	pg	12750	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 SEP 2021