



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 12

Catalog Number: C0092 -2

Lot Number: A00U0278

Expiration Date: 30 SEP 2027

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	11900	pg	47600	pg/mL
IL-17A/F	13625	pg	54500	pg/mL
IP-10	5350	pg	21400	pg/mL
MCP-1	1613	pg	6450	pg/mL
MIP-1 α	2538	pg	10150	pg/mL
MIP-1 β	34125	pg	136500	pg/mL
MIP-2	3038	pg	12150	pg/mL
MIP-3 α	2925	pg	11700	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		26 OCT 2022