



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 12

Catalog Number: C0092 -2

Lot Number: A00U0263

Expiration Date: 31 MAR 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	12500	pg	50000	pg/mL
IL-17A/F	13625	pg	54500	pg/mL
IP-10	6338	pg	25350	pg/mL
MCP-1	1363	pg	5450	pg/mL
MIP-1 α	1950	pg	7800	pg/mL
MIP-1 β	35125	pg	140500	pg/mL
MIP-2	3200	pg	12800	pg/mL
MIP-3 α	3438	pg	13750	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		02 JUN 2022