



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 12  
**Catalog Number:** C0092-2  
**Lot Number:** A00U0406  
**Expiration Date:** 31 MAR 2030  
**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	9512.5	pg	38050	pg/mL
IL-17A/F	12875	pg	51500	pg/mL
IP-10	6712.5	pg	26850	pg/mL
MCP-1	1625	pg	6500	pg/mL
MIP-1 $\alpha$	2587.5	pg	10350	pg/mL
MIP-1 $\beta$	34625	pg	138500	pg/mL
MIP-2	2962.5	pg	11850	pg/mL
MIP-3 $\alpha$	3062.5	pg	12250	pg/mL

**Reconstitution Volume:** 250  $\mu$ 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
<b>Review/Approval</b>	Karen Hamilla	Quality		21 APR 2025