



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 12

Catalog Number: C0092-2

Lot Number: A00U0129

Expiration Date: 31 OCT 2024

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	9463	pg	37850	pg/mL
IL-17A/F	12400	pg	49600	pg/mL
IP-10	5675	pg	22700	pg/mL
MCP-1	1813	pg	7250	pg/mL
MIP-1 α	2163	pg	8650	pg/mL
MIP-1 β	33000	pg	132000	pg/mL
MIP-2	2275	pg	9100	pg/mL
MIP-3 α	3363	pg	13450	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		22 OCT 2019