



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 16

**Catalog Number:** C0295-2

**Lot Number:** A00U0187

**Expiration Date:** 31 AUG 2023

**Storage Temperature:**  $\leq -70^{\circ}\text{C}$

**Intended Assay Species:** Mouse

### Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Calibrator	Unit of Measure
BAFF	4175	pg	16700	pg/mL
6CKine/CCL21	2338	pg	9350	pg/mL
CD40	9500	pg	38000	pg/mL
BCA-1/BLC	26875	pg	107500	pg/mL
IFN- $\beta$	7750	pg	31000	pg/mL
MDC	5950	pg	23800	pg/mL
MCP-5/CCL12	1209	pg	4835	pg/mL

**Volume:** 250  $\mu\text{L}$

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 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		12 APR 2021