



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 16

**Catalog Number:** C0295-2

**Lot Number:** A00U0120

**Expiration Date:** 31 DEC 2021

**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	4 488	pg	17 950	pg/mL
6CKine/CCL21	2 375	pg	9 500	pg/mL
CD40	9 925	pg	39 700	pg/mL
BCA-1/BLC	26 625	pg	106 500	pg/mL
IFN- $\beta$	7 763	pg	31 050	pg/mL
MDC	6 713	pg	26 850	pg/mL
MCP-5/CCL12	1 215	pg	4 860	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 AUG 2019