



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 16

Catalog Number: C0295-2

Lot Number: A00U0224

Expiration Date: 31 DEC 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	4263	pg	17050	pg/mL
6CKine/CCL21	2075	pg	8300	pg/mL
CD40	10200	pg	40800	pg/mL
BCA-1/BLC	23750	pg	95000	pg/mL
IFN- β	8013	pg	32050	pg/mL
MDC	4813	pg	19250	pg/mL
MCP-5/CCL12	1215	pg	4860	pg/mL

Volume: 250 μL

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
Review/Approval	Karen Hamilla	Quality		14 SEP 2021