



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 16

Catalog Number: C0295-2

Lot Number: A00U0267

Expiration Date: 30 NOV 2024

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	4600	pg	18400	pg/mL
6CKine/CCL21	2275	pg	9100	pg/mL
CD40	10575	pg	42300	pg/mL
BCA-1/BLC	26250	pg	105000	pg/mL
IFN- β	7413	pg	29650	pg/mL
MDC	5213	pg	20850	pg/mL
MCP-5/CCL12	1275	pg	5100	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		29 JUN 2022