



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 21

Catalog Number: C0329-2

Lot Number: A00U0156

Expiration Date: 30 APR 2025

Storage Temperature: 2-8°C

Intended Assay Species: Human

Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
BAFF-R/TNFRSF13C	33250	pg	133000	pg/mL
CD40L (soluble)	3750	pg	15000	pg/mL
FGF (basic)	3025	pg	12100	pg/mL
GITR/TNFRSF18	3200	pg	12800	pg/mL
PD-L1 (epitope 1)	2525	pg	10100	pg/mL
PD-L1 (epitope 2)	2525	pg	10100	pg/mL
PD-L2	86250	pg	345000	pg/mL
PIGF	2215	pg	8860	pg/mL
RANKL/TNFSF11	9100	pg	36400	pg/mL
VEGF-D	3525	pg	14100	pg/mL

Reconstitution 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 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		23 NOV 2020