



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 21
Catalog Number: C0329-2
Lot Number: A00U0457
Expiration Date: 31 AUG 2030
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	35750	pg	143000	pg/mL
CD40L (soluble)	4425	pg	17700	pg/mL
FGF (basic)	3475	pg	13900	pg/mL
GITR/TNFRSF18	3450	pg	13800	pg/mL
PD-L1 (epitope 1)	2925	pg	11700	pg/mL
PD-L1 (epitope 2)	2925	pg	11700	pg/mL
PD-L2	104750	pg	419000	pg/mL
PlGF	2377.5	pg	9510	pg/mL
RANKL/TNFSF11	9925	pg	39700	pg/mL
VEGF-D	3275	pg	13100	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		30 SEP 2025