



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 21
Catalog Number: C0329-2
Lot Number: A00U0303
Expiration Date: 31 JAN 2028
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
BAF-R/TNFRSF13C	34500	pg	138000	pg/mL
CD40L (soluble)	3925	pg	15700	pg/mL
FGF (basic)	2285	pg	9140	pg/mL
GITR/TNFRSF18	3425	pg	13700	pg/mL
PD-L1(epitope 1)	1865	pg	7460	pg/mL
PD-L1(epitope 2)	1865	pg	7460	pg/mL
PD-L2	86750	pg	347000	pg/mL
PIGF	2290	pg	9160	pg/mL
RANKL/TNFSF11	9775	pg	39100	pg/mL
VEGF-D	3325	pg	13300	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 was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		07 MAR 2023