



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 31
Catalog Number: C0718-2
Lot Number: A00U0428
Expiration Date: 31 JAN 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
4-1BBL/TNFSF9	35500	pg	142000	pg/mL
CD40/TNFRSF5	3800	pg	15200	pg/mL
FasL	3825	pg	15300	pg/mL
HAVCR1/KIM-1	7825	pg	31300	pg/mL
TACI/TNFRSF13B	159250	pg	637000	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		19 MAR 2025