



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 22
Catalog Number: C0330-2
Lot Number: A00U0480
Expiration Date: 30 JUN 2031
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
BCMA/TNFRSF17	1755	pg	7020	pg/mL
GP130 (soluble)	572500	pg	2290000	pg/mL
HAVCR2/TIM-3	26250	pg	105000	pg/mL
Tie-2	73250	pg	293000	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		28 JUL 2026