



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 22
Catalog Number: C0330-2
Lot Number: A00U0352
Expiration Date: 31 OCT 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
BCMA/TNFRSF17	1625	pg	6500	pg/mL
GP130 (soluble)	532500	pg	2130000	pg/mL
HAVCR2/TIM-3	23750	pg	95000	pg/mL
Tie-2	79000	pg	316000	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		15 DEC 2023