



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 22  
**Catalog Number:** C0330-2  
**Lot Number:** A00U0415  
**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
BCMA/TNFRSF17	1647.5	pg	6590	pg/mL
GP130 (soluble)	557500	pg	2230000	pg/mL
HAVCR2/TIM-3	24700	pg	98800	pg/mL
Tie-2	76000	pg	304000	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
<b>Review/Approval</b>	Karen Hamilla	Quality		11 MAR 2025