

**CERTIFICATE OF ANALYSIS**

Product Description: Calibrator 27
Catalog Number: C0410-2
Lot Number: A00U0434
Expiration Date: 31 MAY 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
APRIL/TNFSF13	145750	pg	583000	pg/mL
E-Selectin	487500	pg	1950000	pg/mL
Galectin-9	13050	pg	52200	pg/mL
HVEM/TNFRSF14	12300	pg	49200	pg/mL
MIG	16500	pg	66000	pg/mL
Pentraxin 3	105500	pg	422000	pg/mL
RAGE (soluble)	4725	pg	18900	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		23 JUL 2025