



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 27
Catalog Number: C0410-2
Lot Number: A00U0466
Expiration Date: 28 FEB 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
APRIL/TNFSF13	132500	pg	530000	pg/mL
E-Selectin	637500	pg	2550000	pg/mL
Galectin-9	13675	pg	54700	pg/mL
HVEM/TNFRSF14	12075	pg	48300	pg/mL
MIG	14325	pg	57300	pg/mL
Pentraxin 3	105750	pg	423000	pg/mL
RAGE (soluble)	4775	pg	19100	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		03 APR 2026