



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 27  
**Catalog Number:** C0410-2  
**Lot Number:** A00U0389  
**Expiration Date:** 31 AUG 2029  
**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	129250	pg	517000	pg/mL
E-Selectin	745000	pg	2980000	pg/mL
Galectin-9	13200	pg	52800	pg/mL
HVEM/TNFRSF14	20050	pg	80200	pg/mL
MIG	15025	pg	60100	pg/mL
Pentraxin 3	163750	pg	655000	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
<b>Review/Approval</b>	Karen Hamilla	Quality		01 OCT 2024