



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 27

**Catalog Number:** C0410-2

**Lot Number:** A00U0265

**Expiration Date:** 30 APR 2027

**Storage Temperature:** 2-8°C

**Intended Assay Species:** Human

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
APRIL/TNFSF13	138250	pg	553000	pg/mL
E-Selectin	775000	pg	3100000	pg/mL
Galectin-9	22300	pg	89200	pg/mL
HVEM/TNFRSF14	21900	pg	87600	pg/mL
MIG	12625	pg	50500	pg/mL
Pentraxin 3	177750	pg	711000	pg/mL
RAGE (soluble)	8250	pg	33000	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		03 JUN 2022