



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 27
Catalog Number: C0410-2
Lot Number: A00U0329
Expiration Date: 31 MAY 2028
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	130000	pg	520000	pg/mL
E-Selectin	750000	pg	3000000	pg/mL
Galectin-9	22275	pg	89100	pg/mL
HVEM/TNFRSF14	21775	pg	87100	pg/mL
MIG	13825	pg	55300	pg/mL
Pentraxin 3	174500	pg	698000	pg/mL
RAGE (soluble)	8600	pg	34400	pg/mL

Reconstitution Volume: 250µL 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 was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		29 JUN 2023