



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 24
Catalog Number: C0351-2
Lot Number: A00U0465
Expiration Date: 31 JAN 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
DPPIV	18500	pg	74000	pg/mL
ICAM-1	2122.5	pg	8490	pg/mL
SAA	20575	pg	82300	pg/mL
SHBG	65250	pg	261000	pg/mL
VCAM-1	59000	pg	236000	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