



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 25
Catalog Number: C0352-2
Lot Number: A00U0427
Expiration Date: 31 JAN 2030
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
Carb. Anh. 1	18400	pg	73600	pg/mL
Complement factor D	9625	pg	38500	pg/mL
CRP	17825	pg	71300	pg/mL
Cystatin C	66750	pg	267000	pg/mL
Factor VII	14475	pg	57900	pg/mL
NGAL/LCN2	12350	pg	49400	pg/mL
sTfR-1	9175	pg	36700	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 Hamila	Quality		04 MAR 2025