



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 26
Catalog Number: C0353-2
Lot Number: A00U0393
Expiration Date: 31 OCT 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
A2M	387500	pg	1550000	pg/mL
Adinopectin	12575	pg	50300	pg/mL
ApoA1	62750	pg	251000	pg/mL
ApoC3	13500	pg	54000	pg/mL
Complement C9	55750	pg	223000	pg/mL
RBP4	16375	pg	65500	pg/mL
Serpin A1	109750	pg	439000	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	Michelle Likiam	Quality		02 DEC 2024