



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 14
Catalog Number: C0272-2
Lot Number: A00U0350
Expiration Date: 31 OCT 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
BAFF	1165	pg	4660	pg/mL
β -NGF	1082.5	pg	4330	pg/mL
BDNF	4300	pg	17200	pg/mL
FGF-23	6925	pg	27700	pg/mL
LH	66500	μ IU	266000	μ IU/mL
FSH	172250	μ IU	689000	μ IU/mL

Reconstitution Volume: 250 μ L

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		06 DEC 2023