



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 14
Catalog Number: C0272-2
Lot Number: A00U0386
Expiration Date: 31 AUG 2029
Storage Temperature: 2-8°C
Intended Assay Species:


Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
BAFF	1272.5	pg	5090	pg/mL
β -NGF	1135	pg	4540	pg/mL
BDNF	4600	pg	18400	pg/mL
FGF-23	5825	pg	23300	pg/mL
LH	70500	μ IU	282000	μ IU/mL
FSH	192500	μ IU	770000	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		25 SEP 2024