



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 14
Catalog Number: C0272-2
Lot Number: A00U0407
Expiration Date: 31 MAR 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
BAFF	1210	pg	4840	pg/mL
β-NGF	1030	pg	4120	pg/mL
BDNF	3725	pg	14900	pg/mL
FGF-23	6475	pg	25900	pg/mL
LH	67500	μIU	270000	μIU/mL
FSH	174250	μIU	697000	μIU/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		29 MAY 2025