



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 14
Catalog Number: C0272-2
Lot Number: A00U0166
Expiration Date: 31 AUG 2025
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	1173	pg	4690	pg/mL
β -NGF	1068	pg	4270	pg/mL
BDNF	4350	pg	17400	pg/mL
FGF-23	6050	pg	24200	pg/mL
LH	68000	μ IU	272000	μ IU/mL
FSH	160750	μ IU	643000	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		28 AUG 2020