



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 14

Catalog Number: C0272-2

Lot Number: A00U0218

Expiration Date: 30 APR 2026

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	1190	pg	4760	pg/mL
β -NGF	1113	pg	4450	pg/mL
BDNF	4225	pg	16900	pg/mL
FGF-23	6375	pg	25500	pg/mL
LH	68250	μ IU	273000	μ IU/mL
FSH	168250	μ IU	673000	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		13 SEP 2021