



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 14  
**Catalog Number:** C0272-2  
**Lot Number:** A00U0366  
**Expiration Date:** 30 APR 2029  
**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    | 1277.5                          | pg              | 5110  | pg/mL           |
| β-NGF   | 1230                            | pg              | 4920  | pg/mL           |
| BDNF    | 5200                            | pg              | 20800   | pg/mL           |
| FGF-23  | 6800                            | pg              | 27200   | pg/mL           |
| LH      | 73250                           | μIU             | 293000  | μIU/mL          |
| FSH     | 212250                          | μIU             | 849000  | 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        |
|------------------------|---------------|----------|-----------|-------------|
| <b>Review/Approval</b> | Karen Hamilla | Quality  |           | 05 JUN 2024 |