



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 14

**Catalog Number:** C0272-2

**Lot Number:** A00U0281

**Expiration Date:** 31 AUG 2027

**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	1260	pg	5040	pg/mL
$\beta$ -NGF	1215	pg	4860	pg/mL
BDNF	4925	pg	19700	pg/mL
FGF-23	6575	pg	26300	pg/mL
LH	71000	$\mu$ IU	284000	$\mu$ IU/mL
FSH	176250	$\mu$ IU	705000	$\mu$ IU/mL

**Reconstitution Volume:** 250  $\mu$ 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
<b>Review/Approval</b>	Karen Hamilla	Quality		05 OCT 2022