



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 14

**Catalog Number:** C0272-2

**Lot Number:** A00U0067

**Expiration Date:** 31 JUL 2023

**Storage Condition:** 2-8°C

Analyte Description	Amount of Analyte in Calibrator	Units	Stock Concentration of Reconstituted Calibrator (Vol = 250 µL)	Units
hu BAFF	1 023	pg	4 090	pg/mL
hu β-NGF	1 090	pg	4 360	pg/mL
hu BDNF	4 100	pg	16 400	pg/mL
hu FGF-23	6 325	pg	25 300	pg/mL
hu LH	67 750	µIU	271 000	µIU/mL
hu FSH	162 000	µIU	648 000	µIU/mL

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.

**Statement:** The above calibrator lot is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	30 AUG 2018