



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 13

**Catalog Number:** C0271-2

**Lot Number:** A00U0066

**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 GIP (total)	27 800	pg	111 000	pg/mL
hu GIP (inactive)	27 800	pg	111 000	pg/mL
hu GLP-1 (total)	1.82	pmol	7 270	pM
hu GLP-1 (inactive)	1.82	pmol	7 270	pM
hu FGF-21	18 000	pg	71 800	pg/mL
hu Ghrelin (total)	11 000	pg	44 100	pg/mL
hu Glucagon	0.400	pmol	1 600	pM
hu Leptin	121 000	pg	483 000	pg/mL
hu PP	4 300	pg	17 200	pg/mL
hu PYY (total)	5 700	pg	22 800	pg/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
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	30 AUG 2018