



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 13

**Catalog Number:** C0271-2

**Lot Number:** A00U0122

**Expiration Date:** 31 AUG 2024

**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
GIP (total)	31 500	pg	126 000	pg/mL
GIP (inactive)	31 500	pg	126 000	pg/mL
GLP-1 (total)	1.84	pmol	7 350	pM
GLP-1 (inactive)	1.84	pmol	7 350	pM
FGF-21	20 450	pg	81 800	pg/mL
Ghrelin (total)	12 950	pg	51 800	pg/mL
Glucagon	0.393	pmol	1 570	pM
Leptin	122 000	pg	488 000	pg/mL
PP	4 325	pg	17 300	pg/mL
PYY (total)	5 700	pg	22 800	pg/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		20 SEP 2019