



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 13  
**Catalog Number:** C0271-2  
**Lot Number:** A00U0414  
**Expiration Date:** 28 FEB 2030  
**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)	35000	pg	140000	pg/mL
GLP-1 (total)	1.7975	pmol	7190	pM
FGF-21	27000	pg	108000	pg/mL
Ghrelin (total)	13550	pg	54200	pg/mL
Glucagon	0.395	pmol	1580	pM
Leptin	115500	pg	462000	pg/mL
PP	4250	pg	17000	pg/mL
PYY (total)	6075	pg	24300	pg/mL
GIP (inactive)	35000	pg	140000	pg/mL
GLP-1 (inactive)	1.7975	pmol	7190	pM

**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		10 APR 2025