



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 13
Catalog Number: C0271-2
Lot Number: A00U0454
Expiration Date: 31 AUG 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 (inactive)	27000	pg	108000	pg/mL
GIP (total)	27000	pg	108000	pg/mL
GLP-1 (inactive)	1.76	pmol	7040	pM
GLP-1 (total)	1.76	pmol	7040	pM
FGF-21	19350	pg	77400	pg/mL
Ghrelin (total)	14850	pg	59400	pg/mL
Glucagon	0.375	pmol	1500	pM
Leptin	112250	pg	449000	pg/mL
PP	4300	pg	17200	pg/mL
PYY (total)	6200	pg	24800	pg/mL

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
Review/Approval	Karen Hamilla	Quality		25 SEP 2025