



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 13

Catalog Number: C0271-2

Lot Number: A00U0284

Expiration Date: 31 AUG 2027

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)	34750	pg	139000	pg/mL
GLP-1 (total)	1.95	pmol	7800	pM
FGF-21	23150	pg	92600	pg/mL
Ghrelin (total)	15650	pg	62600	pg/mL
Glucagon	0.390	pmol	1560	pM
Leptin	124250	pg	497000	pg/mL
PP	4425	pg	17700	pg/mL
PYY (total)	5925	pg	23700	pg/mL
GIP (inactive)	34750	pg	139000	pg/mL
GLP-1 (inactive)	1.95	pmol	7800	pM

Reconstitution Volume: 250 μ 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
Review/Approval	Karen Hamilla	Quality		13 OCT 2022