



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 13  
**Catalog Number:** C0271-2  
**Lot Number:** A00U0344  
**Expiration Date:** 30 SEP 2028  
**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)	30250	pg	121000	pg/mL
GLP-1 (total)	2.045	pmol	8180	pM
FGF-21	20400	pg	81600	pg/mL
Ghrelin (total)	15700	pg	62800	pg/mL
Glucagon	0.3875	pmol	1550	pM
Leptin	121500	pg	486000	pg/mL
PP	4425	pg	17700	pg/mL
PYY(Total)	5850	pg	23400	pg/mL
GLP-1(inactive)	2.045	pmol	8180	pM
GIP (inactive)	30250	pg	121000	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 was manufactured and tested according to MSD documents. This lot meets MSD's specifications. 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		02 NOV 2023