



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 18  
**Catalog Number:** C0292-2  
**Lot Number:** A00U0306  
**Expiration Date:** 31 JAN 2028  
**Storage Temperature:** 2-8°C  
**Intended Assay Species:** Mouse/Rat

**Concentration Assignment:**

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
GLP-1 (total)	2.069	pmol	8275	pM
GLP-1 (inactive)	2.069	pmol	8275	pM
FGF-21	33500	pg	134000	pg/mL
BDNF	5713	pg	22850	pg/mL
Ghrelin (total)	23838	pg	95350	pg/mL
PYY (total)	9200	pg	36800	pg/mL
C-peptide	288750	pg	1155000	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		29 SEP 2023