



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 18  
**Catalog Number:** C0292-2  
**Lot Number:** A00U0324  
**Expiration Date:** 31 JUL 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)	2.23	pmol	8920	pM
FGF-21	24150	pg	96600	pg/mL
BDNF	6175	pg	24700	pg/mL
Ghrelin (total)	21150	pg	84600	pg/mL
PYY (total)	10000	pg	40000	pg/mL
C-Peptide	292500	pg	1170000	pg/mL
GLP-1 (inactive)	2.23	pmol	8920	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 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		22 SEP 2023