



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 18
Catalog Number: C0292-2
Lot Number: A00U0416
Expiration Date: 28 FEB 2030
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)	2.0125	pmol	8050	pM
GLP-1 (Inactive)	2.0125	pmol	8050	pM
FGF-21	34500	pg	138000	pg/mL
BDNF	6775	pg	27100	pg/mL
Ghrelin (total)	24725	pg	98900	pg/mL
PYY (total)	9950	pg	39800	pg/mL
C-peptide	305000	pg	1220000	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		24 APR 2025