



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 18
Catalog Number: C0292-2
Lot Number: A00U0374
Expiration Date: 30 APR 2029
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)	2.3775	pmol	9510	pM
GLP-1 (inactive)	2.3775	pmol	9510	pM
FGF-21	35250	pg	141000	pg/mL
BDNF	6700	pg	26800	pg/mL
Ghrelin (total)	25000	pg	100000	pg/mL
PYY (total)	9500	pg	38000	pg/mL
C-peptide	300000	pg	1200000	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		28 MAY 2024