



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 18
Catalog Number: C0292-2
Lot Number: A00U0121
Expiration Date: 31 JUL 2024
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)	2.255	pmol	9 020	pM
GLP-1 (inactive)	2.255	pmol	9 020	pM
FGF-21	29 000	pg	116 000	pg/mL
BDNF	5 250	pg	21 000	pg/mL
Ghrelin (total)	25 250	pg	101 000	pg/mL
C-Peptide	322 500	pg	1 290 000	pg/mL
PYY(total)	10 200	pg	40 800	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		12 AUG 2019