



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 18
Catalog Number: C0292-2
Lot Number: A00U0247
Expiration Date: 31 MAR 2027
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)	2.230	pmol	8920	pM
GLP-1 (inactive)	2.230	pmol	8920	pM
FGF-21	29750	pg	119000	pg/mL
BDNF	5850	pg	23400	pg/mL
Ghrelin (total)	26250	pg	105000	pg/mL
PYY (total)	10250	pg	41000	pg/mL
C-peptide	307500	pg	1230000	pg/mL

Reconstitution Volume: 250 μ L

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		26 APR 2022