



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 18
Catalog Number: C0292-2
Lot Number: A00U0195
Expiration Date: 31 JAN 2026
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
BDNF	5525	pg	22100	pg/mL
FGF-21	29250	pg	117000	pg/mL
Ghrelin (total)	22575	pg	90300	pg/mL
GLP-1 (total)	2	pmol	8880	pM
GLP-1 (inactive)	2	pmol	8880	pM
C-peptide	325000	pg	1300000	pg/mL
PYY (total)	10050	pg	40200	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		14 MAY 2021