



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 19
Catalog Number: C0293-2
Lot Number: A00U0186
Expiration Date: 30 NOV 2025
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
Insulin	15900	μIU	63600	μIU/mL
Leptin	141750	pg	567000	pg/mL
Glucagon	398	pmol	1590	pM

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		20 JAN 2021