



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 19
Catalog Number: C0293-2
Lot Number: A00U0321
Expiration Date: 31 MAR 2028
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
Glucagon	0.3475	pmol	1390	pM
Insulin	14775	μIU	59100	μIU/mL
Leptin	140750	pg	563000	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 was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		09 MAY 2023