

## Page 1 of 1

**Product Description:** Ghrelin (active) Calibrator

Catalog Number:C016K-2Lot Number:A00U0282Expiration Date31 JUL 2027

**Storage Temperature:** 2-8°C

**Intended Assay Species:** Human, Mouse Rat

**Concentration Assignment:** 

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Ghrelin (active)	39750	pg	159000	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	Laur Hamille	30 OCT 2024

COA-16938-C QA-FM-184-J