



## CERTIFICATE OF ANALYSIS

COA – 05028

**Product Description:** MSD<sup>®</sup> Glucagon Calibrator

**Catalog #/Size:** C00HC-2 / 1 Plate

**Lot Number:** A0010347

**Expiration Date:** 31 OCT 2016

**Statement:** This product was manufactured and tested according to approved MSD documents. This lot meets MSD's specifications.

Volume: 15 µL

Concentration: 1 µg/mL

Storage buffer: NA

Storage conditions: ≤-70° C

**The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.**

	<b>Name</b>	<b>Function</b>	<b>Signature</b>	<b>Date</b>
<b>Preparer</b>	Rolanda Spratley	Biologics/Chemistry		11 DEC 2014
<b>Review/Approval</b>	Anastasia Belilovets	Quality Control		11 DEC 2014