



**CERTIFICATE OF ANALYSIS**

Page 1 of 1

**Product Description:** GLP-1 (7-36)amide Calibrator

**Catalog Number:** C00FA-2

**Lot Number:** A0010446

**Expiration Date:** 30 SEP 2020

**Storage Temperature:**  $\leq -70^{\circ}\text{C}$

**Concentration of Materials Tested:**

	Assay
	<b>GLP-1 (7-36)amide</b>
Calibrator $\mu\text{g}/\text{mL}$	1.00

Volume: 15  $\mu\text{L}$

**Statement:** This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	17 DEC 2018