



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: GLP-1 (active) Calibrator

Catalog Number: C0266-1

Lot Number: A0080155

Expiration Date: 30 NOV 2022

Storage Temperature: 2-8°C

Concentration of Materials Tested:

	Assay
	GLP-1 (active)
Reconstituted Calibrator (pM)	276

Reconstitution Volume: 1000 µL

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	15 NOV 2017