



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: GLP-1 (active) Calibrator
Catalog Number: C0266-1
Lot Number: A0080195
Expiration Date: 31 JAN 2024
Storage Temperature: 2-8°C

Concentration of Materials Tested:

	Assay
	GLP-1 (active)
Reconstitution Calibrator (pM)	267

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	Kayla Turner	Quality Assurance		30 Jan 2019