



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: GLP-1 (total) Calibrator
Catalog Number: C003P-2
Lot Number: A0080194
Expiration Date: 31 DEC 2023
Storage Temperature: 2-8°C

Concentration of Materials Tested:

	Assay
	GLP-1 (total)
Reconstitution Calibrator (pM)	195

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>	29 JAN 2019