



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: GLP-1 (total) Calibrator

Catalog Number: C003P-2

Lot Number: A0080160

Expiration Date: 31 JAN 2023

Storage Temperature: 2-8°C

Concentration of Materials Tested:

	Assay
	GLP-1 (total)
Calibrator (pM)	194

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>	23 MAR 2018