



CERTIFICATE OF ANALYSIS

COA - 07994

Product Description: MSD® Human Sclerostin Kit

Catalog #/Size: K151HGC-Series

Lot Number: K0050735

Expiration Date: N/A (See expiration date on specific components below)

Statement: This product was manufactured and tested according to approved MSD documents. This lot meets MSD's specifications.

Kit Contents:

Plate Lot Number: Z0055570

Plate Lot Catalog Number: N75146A

Plate Lot Expiration Date: 31 AUG 2018

Antibody blend Lot Number: D0050669

Antibody blend Catalog Number: D21HG

Antibody blend Expiration Date: 31 JUL 2018

Calibrator blend Lot Number: A0020189

Calibrator blend Catalog Number: C0146

Calibrator blend Expiration Date: 30 SEP 2018

Specifications:

Release method: Coating uniformity

Mean Intraplate CV: $\leq 15\%$

Mean Interplate CV: $\leq 18\%$

No more than one plate with CV $> 18\%$

No plate with CV $> 20\%$

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Preparer	Xiomara Talavera	Quality Control		14 MAR 2017
Review/Approval	Vanessa Castleberry	Quality Assurance		15 MAR 2017