



**CERTIFICATE OF ANALYSIS**

Page 1 of 1

**Product Description:** MSD® Human Sclerostin Kit

**Catalog Number:** K151HGC - Series

**Lot Number:** K0050830

**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					
<b>Lot #</b>	Z0056301	<b>Catalog #</b>	N75146A	<b>Expiration Date:</b>	31 DEC 2020

Antibody Blend	
<b>Lot Number:</b>	D0050693
<b>Catalog Number:</b>	D21HG
<b>Expiration Date:</b>	31 MAY 2019
<b>Storage Temperature:</b>	2-8°C

Calibrator Blend	
<b>Lot Number:</b>	A0020247
<b>Catalog Number:</b>	C0146
<b>Expiration Date:</b>	31 DEC 2020
<b>Storage Temperature:</b>	≤ -70°C

Specifications:

Release method:

Coating Uniformity	Specifications
Mean Intraplate CV	≤ 15%
Mean Interplate CV	≤ 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
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	28 NOV 2018