



## CERTIFICATE OF ANALYSIS

COA - 04811

**Product Description:** MSD<sup>®</sup> Human Osteopontin Kit

**Catalog #/Size:** K151HJC-1 / 1 Plate; K151HJC-2 / 5 Plates; K151HJC-3 / 20 Plates

**Lot Number:** K0050677

**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: Z0055046

Plate Lot Catalog Number: N75147A

Plate Lot Expiration Date: 31 JAN 2017

Antibody blend Lot Number: D0050604

Antibody blend Catalog Number: D21HJ

Antibody blend Expiration Date: 31 DEC 2016

Calibrator blend Lot Number: A0020162

Calibrator blend Catalog Number: C0147

Calibrator blend Expiration Date: 30 NOV 2016

### 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
<b>Preparer</b>	Rolanda Spratley	Biologics/Chemistry		10 FEB 2015
<b>Review/Approval</b>	Anastasia Belilovets	Quality Assurance		12 FEB 2015