



## CERTIFICATE OF ANALYSIS

COA - 07798

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

**Catalog #/Size:** K150JLD - Series

**Lot Number:** K00E0312

**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: Z00E0247

Plate Lot Catalog Number: N45203A

Plate Lot Expiration Date: 30 APR 2019

Antibody blend Lot Number: D00E0290

Antibody blend Catalog Number: D20JL

Antibody blend Expiration Date: 30 APR 2019

Calibrator blend Lot Number: A00E0154

Calibrator blend Catalog Number: C0203

Calibrator blend Expiration Date: 31 MAY 2019

### 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.**

	<b>Name</b>	<b>Function</b>	<b>Signature</b>	<b>Date</b>
<b>Preparer</b>	Xiomara Talavera	Quality Control		25 JAN 2017
<b>Review/Approval</b>	Vanessa Castleberry	Quality Assurance		27 JAN 2017