



CERTIFICATE OF ANALYSIS

COA -07072

**Product Description:** MSD® Human/NHP IgG Kit

**Catalog #/Size:** K150JLD-1 / 1 Plate; K150JLD-2 / 5 Plates; K150JLD-4 / 25 Plates

**Lot Number:** K00E0266

**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: Z00E0213
- Plate Lot Catalog Number: N45203A
- Plate Lot Expiration Date: 31 AUG 2018
- Antibody blend Lot Number: D00E0245
- Antibody blend Catalog Number: D20JL
- Antibody blend Expiration Date: 31 AUG 2018
- Calibrator blend Lot Number: A00E0135
- Calibrator blend Catalog Number: C0203
- Calibrator blend Expiration Date: 31 OCT 2018

Specifications:

- Release method: Coating uniformity
- 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>Preparer</b>	Rolanda Spratley	Biologics/Chemistry		30 JUN 2016
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance		30 JUN 2016