



## CERTIFICATE OF ANALYSIS

COA - 06931

**Product Description:** MSD<sup>®</sup> Isotyping Panel 1 (Human/NHP) Kit

**Catalog #/Size:** K15203D-1 / 1 Plate; K15203D-2 / 5 Plates; K15203D-4 / 25 Plates

**Lot Number:** K00E0251

**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: Z00E0206

Plate Lot Catalog Number: N45203A

Plate Lot Expiration Date: 31 MAY 2018

Antibody blend Lot Number: D00E0234

Antibody blend Catalog Number: D20JJ

Antibody blend Expiration Date: 31 MAY 2018

Antibody blend Lot Number: D00E0235

Antibody blend Catalog Number: D20JL

Antibody blend Expiration Date: 30 JUN 2018

Antibody blend Lot Number: D00E0236

Antibody blend Catalog Number: D20JP

Antibody blend Expiration Date: 31 MAY 2018

Calibrator blend Lot Number: A00E0124

Calibrator blend Catalog Number: C0203

Calibrator blend Expiration Date: 30 APR 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	Rolanda Spratley	Biologics/Chemistry		23 MAY 2016
Review/Approval	Dawn Stollar	Quality Assurance		23 MAY 2016