



CERTIFICATE OF ANALYSIS

COA - 06914

Product Description: MSD[®] Human MIP-4 Kit

Catalog #/Size: K151RLD-1 / 1 Plate; K151RLD-2 / 5 Plates; K151RLD-4 / 25 Plates

Lot Number: K00E0219

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: Z00E0184

Plate Lot Catalog Number: N451RLA

Plate Lot Expiration Date: 30 NOV 2017

Antibody blend Lot Number: D00E0207

Antibody blend Catalog Number: D21RL

Antibody blend Expiration Date: 30 NOV 2017

Calibrator blend Lot Number: A00E0112

Calibrator blend Catalog Number: C01RL

Calibrator blend Expiration Date: 30 NOV 2017

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		18 MAY 2016
Review/Approval	Dawn Stollar	Quality Assurance		18 MAY 2016