



CERTIFICATE OF ANALYSIS

COA - 07098

Product Description: MSD[®] Human KIM-1 Kit

Catalog #/Size: K151JHD-1 / 1 Plate; K151JHD-2 / 5 Plates; K151JHD-4 / 25 Plates

Lot Number: K00E0202

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: Z00E0178
- Plate Lot Catalog Number: N451JHA
- Plate Lot Expiration Date: 30 SEP 2017
- Antibody blend Lot Number: D00E0195
- Antibody blend Catalog Number: D21JH
- Antibody blend Expiration Date: 30 SEP 2017
- Calibrator blend Lot Number: A00E0104
- Calibrator blend Catalog Number: C01JH
- Calibrator blend Expiration Date: 30 SEP 2017

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
Preparer	Rolanda Spratley	Biologics/Chemistry		11 JUL 2016
Review/Approval	Vanessa Castleberry	Quality Assurance		11 JUL 2016