



CERTIFICATE OF ANALYSIS

COA - 07380

Product Description: MSD[®] Human IL-6R Ultra-Sensitive Kit

Catalog #/Size: K151ALC-1 / 1 Plate; K151ALC-2 / 5 Plates; K151ALC-4 / 25 Plates

Lot Number: K0035124

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: Z0046053

Plate Lot Catalog Number: L451ALA

Plate Lot Expiration Date: 31 AUG 2018

Antibody blend Lot Number: D0033600

Antibody blend Catalog Number: D21AL

Antibody blend Expiration Date: 31 JUL 2018

Antibody source (host/species): Goat

Calibrator blend Lot Number: A0032221

Calibrator blend Catalog Number: C01AL

Calibrator blend Expiration Date: 31 AUG 2018

Calibrator source (host/species): Sf 21

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	Xiomara Talavera	Quality Control	<i>X Talavera</i>	06 OCT 2016
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	07 OCT 2016