



CERTIFICATE OF ANALYSIS

COA - 04732

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

Catalog #/Size: K151AKC-1 / 1 Plate; K151AKC-2 / 5 Plates; K151AKC-4 / 25 Plates

Lot Number: K0034133

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: Z0044762

Plate Lot Catalog Number: L451AKA

Plate Lot Expiration Date: 30 NOV 2015

Antibody blend Lot Number: D0032814

Antibody blend Catalog Number: D21AK

Antibody blend Expiration Date: 31 DEC 2015

Antibody source (host/species): Goat

Calibrator blend Lot Number: A0031799

Calibrator blend Catalog Number: C01AK

Calibrator blend Expiration Date: 15 Months From Date of Receipt

Calibrator source (host/species): *E. coli*

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		08 SEP 2014
Review/Approval	Anastasia Belilovets	Quality Control		11 SEP 2014