



CERTIFICATE OF ANALYSIS

COA - 03964

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: K0034046

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: Z0044650

Plate Lot Catalog Number: L451AKA

Plate Lot Expiration Date: 30 JUN 2014

Antibody blend Lot Number: D0032742

Antibody blend Catalog Number: D21AK

Antibody blend Expiration Date: 30 APR 2015

Antibody source (host/species): Rabbit

Calibrator blend Lot Number: A0031762

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: ≤ 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		25 Nov 2013
Review/Approval	Dana Short	Quality Control		26 Nov 2013