



CERTIFICATE OF ANALYSIS

COA - 06663

Product Description: MSD[®] Human ProInflammatory-4 I Tissue Culture Kit

Catalog #/Size: K15009B-1 / 1 Plate; K15009B-2 / 5 Plates; K15009B-4 / 25 Plates

Lot Number: K0034826

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: Z0045578
- Plate Lot Catalog Number: N45009B
- Plate Lot Expiration Date: 30 JUN 2017
- Antibody blend Lot Number: D0033346
- Antibody blend Catalog Number: D2009
- Antibody blend Expiration Date: 30 JUN 2017
- Antibody source (host/species): IL-1 β , IL-6, TNF- α : Goat
IFN- γ : Mouse
- Calibrator blend Lot Number: A0032116
- Calibrator blend Catalog Number: C01XX
- Calibrator blend Expiration Date: 30 APR 2017
- 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		22 MAR 2016
Review/Approval	Dawn Stollar	Quality Assurance		22 MAR 2016