



CERTIFICATE OF ANALYSIS

COA - 05037

Product Description: MSD[®] Mouse ProInflammatory 7-Plex Tissue Culture Kit
Catalog #/Size: K15012B-1 / 1 Plate; K15012B-2 / 5 Plates; K15012B-4 / 25 Plates
Lot Number: K0034300

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: Z0044965
Plate Lot Catalog Number: N75012B
Plate Lot Expiration Date: 29 FEB 2016
Antibody blend Lot Number: D0032958
Antibody blend Catalog Number: D2012
Antibody blend Expiration Date: 31 MAR 2016
Antibody source (host/species): mIL-1 β , mIL-12p70, mIL-6, mIL-10, mKC: Goat
mIFN γ , mTNF α : Rat
Calibrator blend Lot Number: A0031879
Calibrator blend Catalog Number: C0012
Calibrator blend Expiration Date: 31 MAR 2016
Calibrator source (host/species): mIL-1 β , mIL-6, mIL-10, mKC, mIFN γ , mTNF α : *E. coli*
mIL-12p70: *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	Rolanda Spratley	Biologics/Chemistry		16 DEC 2014
Review/Approval	Anastasia Belilovets	Quality Control		18 DEC 2014