



CERTIFICATE OF ANALYSIS

COA - 06273

Product Description: MSD[®] Human TH1/TH2 7-Plex Tissue Culture Kit

Catalog #/Size: K15011B-1 / 1 Plate; K15011B-2 / 5 Plates; K15011B-4 / 25 Plates

Lot Number: K0034872

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: Z0045645

Plate Lot Catalog Number: N75011B

Plate Lot Expiration Date: 31 AUG 2017

Antibody blend Lot Number: D0033386

Antibody blend Catalog Number: D2011

Antibody blend Expiration Date: 31 AUG 2017

Antibody source (host/species): hIFN γ , hIL-2, hIL-5, hIL-12p40/70: Mouse

hIL-13: Goat

hIL-4, hIL-10: Rat

Calibrator blend Lot Number: A0032134

Calibrator blend Catalog Number: C01XX

Calibrator blend Expiration Date: 31 OCT 2017

Calibrator source (host/species): hIFN γ , hIL-2, hIL-4, hIL-13: *E. coli*

hIL-5, hIL-10, hIL-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		08 DEC 2015
Review/Approval	Dawn Stollar	Quality Assurance		08 DEC 2015