



CERTIFICATE OF ANALYSIS

COA - 07367

Kit Description: MSD® Human TH1/TH2 10-Plex Tissue Culture Kit
Catalog #/Size: K15010B-1 / 1 Plate; K15010B-2 / 5 Plates; K15010B-4 / 25 Plates, K15010B-5 / 250 Plates
Lot Number: K0034517
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: Z0045103
Plate Lot Catalog Number: N05010B
Plate Lot Expiration Date: 30 JUN 2016
Antibody blend Lot Number: D0033033
Antibody blend Catalog Number: D2010
Antibody blend Expiration Date: 31 JUL 2016
Antibody source (host/species): hIFNγ, hIL-2, hIL-4, hIL-12p70: Mouse
hIL-1β, hIL-8, hIL-13, hTNFα: Goat
hIL-5, hIL-10: Rat
Calibrator blend Lot Number: A0031942
Calibrator blend Catalog Number: C0010
Calibrator blend Expiration Date: 31 AUG 2016
Calibrator source (host/species): hIL-1β/ hIL-1F2, hIL-2, hIL-4, hIL-8, hIL-13,
hIFNγ, hNFα: E. coli
hIL-5, hIL-12p70, hIL-10: Sf 21

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.

Table with 5 columns: Name, Function, Signature, Date. Rows include Preparer (Xiomara Talavera) and Review/Approval (Dawn Stollar).