



CERTIFICATE OF ANALYSIS

COA - 07037

Kit Description: MSD® Mouse TH1/TH2 9-Plex Tissue Culture Kit
Catalog #/Size: K15013B-1 / 1 Plate; K15013B-2 / 5 Plates; K15013B-4 / 25 Plates
Lot Number: K0034977
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: Z0045813
Plate Lot Catalog Number: N05013B
Plate Lot Expiration Date: 31 JAN 2018
Antibody blend Lot Number: D0033475
Antibody blend Catalog Number: D2013
Antibody blend Expiration Date: 31 DEC 2017
Antibody source (host/species): mIFNγ, mIL-2, mIL-4, mIL-5, mTNFα: Rat
mIL-1β, mKC/GRO, mIL-10, mIL-12total: Goat
Calibrator blend Lot Number: A0032155
Calibrator blend Catalog Number: C0013
Calibrator blend Expiration Date: 31 DEC 2017
Calibrator source (host/species): mIL-1β, mIL-2, mIL-4, mIL-10, mKC/GRO, mIFNγ, mTNFα: E. coli
mIL-5, mIL-12total: 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 (Rolanda Spratley) and Review/Approval (Vanessa Castleberry).