



CERTIFICATE OF ANALYSIS

COA - 06280

Product Description: MSD[®] Mouse/Rat EPO Serum/Plasma Kit
Catalog #/Size: K150BQC-1 / 1 Plate; K150BQC-2 / 5 Plates; K150BQC-3 / 20 Plates
Lot Number: K0013658
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: Z0055091
Plate Lot Catalog Number: N45123B
Plate Lot Expiration Date: 31 DEC 2016
Antibody blend Lot Number: D0013235
Antibody blend Catalog Number: D20BQ
Antibody blend Expiration Date: 31 DEC 2016
Calibrator blend Lot Number: A0010354
Calibrator blend Catalog Number: C02BQ
Calibrator blend Expiration Date: 31 OCT 2016
Calibrator blend Lot Number: A0010356
Calibrator blend Catalog Number: C03BQ
Calibrator blend Expiration Date: 05 NOV 2016

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.

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biologics/Chemistry		04 DEC 2015
Review/Approval	Dawn Stollar	Quality Assurance		09 DEC 2015