



CERTIFICATE OF ANALYSIS

COA - 07250

Product Description: MSD® Isotyping Panel 1 (Human/NHP) Kit
Catalog #/Size: K15203D-1 / 1 Plate; K15203D-2 / 5 Plates; K15203D-4 / 25 Plates
Lot Number: K00E0264
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: Z00E0213
 Plate Lot Catalog Number: N45203A
 Plate Lot Expiration Date: 31 AUG 2018
 Antibody blend Lot Number: D00E0244
 Antibody blend Catalog Number: D20JJ
 Antibody blend Expiration Date: 31 AUG 2018
 Antibody blend Lot Number: D00E0245
 Antibody blend Catalog Number: D20JL
 Antibody blend Expiration Date: 31 AUG 2018
 Antibody blend Lot Number: D00E0246
 Antibody blend Catalog Number: D20JP
 Antibody blend Expiration Date: 31 AUG 2018
 Calibrator blend Lot Number: A00E0135
 Calibrator blend Catalog Number: C0203
 Calibrator blend Expiration Date: 31 OCT 2018

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	Xiomara Talavera	Quality Control	<i>Xiomara Talavera</i>	26 AUG 2016
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	29 AUG 2016