



CERTIFICATE OF ANALYSIS

COA - 07865

Product Description: MSD® Isotyping Panel 1 (Human/NHP) Kit

Catalog #/Size: K15203D

Lot Number: K00E0290

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: Z00E0234
- Plate Lot Catalog Number: N45203A
- Plate Lot Expiration Date: 31 DEC 2018
- Antibody blend Lot Number: D00E0270
- Antibody blend Catalog Number: D20JJ
- Antibody blend Expiration Date: 31 DEC 2018
- Antibody blend Lot Number: D00E0271
- Antibody blend Catalog Number: D20JL
- Antibody blend Expiration Date: 31 DEC 2018
- Antibody blend Lot Number: D00E0272
- Antibody blend Catalog Number: D20JP
- Antibody blend Expiration Date: 31 DEC 2018
- Calibrator blend Lot Number: A00E0147
- Calibrator blend Catalog Number: C0203
- Calibrator blend Expiration Date: 31 DEC 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>	09 FEB 2017
Review/Approval	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	09 FEB 2017