



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: K15203D - Series

Lot Number: K00E0355

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 # | Z00E0278 | Catalog # | N45203A | Expiration Date: | 30 NOV 2019 |

| Antibody Blend | |
|-----------------------------|-------------|
| LotNumber: | D00E0331 |
| Catalog Number: | D20JJ |
| Expiration Date: | 31 JAN 2020 |
| Storage Temperature: | 2-8°C |

| Calibrator Blend | |
|-----------------------------|-------------|
| Lot Number: | A00E0178 |
| Catalog Number: | C0203 |
| Expiration Date: | 31 DEC 2019 |
| Storage Temperature: | ≤ -70°C |

| Antibody Blend | |
|-----------------------------|-------------|
| LotNumber: | D00E0332 |
| Catalog Number: | D20JL |
| Expiration Date: | 31 DEC 2019 |
| Storage Temperature: | 2-8°C |

| Antibody Blend | |
|-----------------------------|-------------|
| LotNumber: | D00E0333 |
| Catalog Number: | D20JP |
| Expiration Date: | 31 DEC 2019 |
| Storage Temperature: | 2-8°C |

Specifications:

Release method:

| Coating Uniformity | Specifications |
|--------------------------------|----------------|
| 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 |
|------------------------|--------------|-------------------|---------------------|-------------|
| Review/Approval | Dawn Stollar | Quality Assurance | <i>Dawn Stollar</i> | 18 OCT 2017 |