



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: sAPPalpha/sAPPbeta Kit

Catalog Number: K15120E - Series

Lot Number: K0014015

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 # | Z0055771 | Catalog # | N45120B | Expiration Date: | 30 SEP 2019 |

| Antibody Blend | |
|----------------------|-------------|
| Lot Number: | D0013622 |
| Catalog Number: | D21EA |
| Expiration Date: | 30 APR 2019 |
| Storage Temperature: | 2-8°C |

| Calibrator Blend | |
|----------------------|-------------|
| Lot Number: | A0010391 |
| Catalog Number: | C01BS |
| Expiration Date: | 31 OCT 2018 |
| Storage Temperature: | ≤ -70°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> | 09 OCT 2017 |