



RFE MEMO

Page 1 of 1

Product Description(s): V-PLEX® and V-PLEX Plus Human TARC Kit

Catalog Number(s) / Size(s): K151NT-Series

Lot Number: K0080771

Based on additional testing, the original date of expiration for this Kit has been extended by 12 months. The expiration date of K0080771 has been extended to 30 JUN 2019.

The expiration dates for the following kit components have also been extended:

| Description | Lot Number | Original Expiration Date | New Expiration Date |
|-----------------------|------------|--------------------------|---------------------|
| Anti-hu TARC Antibody | D0080693 | 30 JUN 2018 | 30 JUN 2019 |
| Diluent 43 | M0430017 | 31 DEC 2018 | 30 JUN 2019 |
| Diluent 3 | M0030070 | 31 JUL 2018 | 30 JUN 2019 |

| Function | Name | Signature | Date |
|-------------------|--------------|---------------------|-------------|
| Quality Assurance | Dawn Stollar | <i>Dawn Stollar</i> | 14 MAR 2018 |

This memo supersedes all other statements regarding date of expiration for this product.

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human TARC Kit
Kit Catalog Numbers: K151NT-series
Kit Lot Number: K0080771
Expiration Date: 30 Jun 2018

Acceptance Data/ Percent Recovery Specifications:

| Spot | Name | Calibrator | | Chemokine panel 1(human) Control 1 | | Chemokine panel 1(human) Control 2 | | Chemokine panel 1(human) Control 3 | |
|------|---------|------------|-------|------------------------------------|-------|------------------------------------|-------|------------------------------------|-------|
| | | Recovery | CV | Recovery | CV | Recovery | CV | Recovery | CV |
| 1 | hu TARC | 80% - 120% | < 20% | 80% - 120%* | < 20% | 80% - 120%* | < 20% | 80% - 120%* | < 20% |

* See Additional Comments.

| Spot | Name | LLOQ | | | ULOQ | | |
|------|---------|-------|------------|-------|-------|------------|-------|
| | | Value | Recovery | CV | Value | Recovery | CV |
| 1 | hu TARC | 3.32 | 80% - 120% | < 20% | 1120 | 80% - 120% | < 20% |

Stock Concentration of Material Tested :

| Spot | Name | Item | Conc | Units |
|------|---------|------------------------------------|------|-------|
| 1 | hu TARC | Calibrator | 1890 | pg/mL |
| | | Chemokine panel 1(human) Control 1 | 1160 | pg/mL |
| | | Chemokine panel 1(human) Control 2 | 188 | pg/mL |
| | | Chemokine panel 1(human) Control 3 | 34.3 | pg/mL |

Controls tested may not be the same lot as shipped.

Standard Curve Data:

| hu TARC - Spot 1 | | |
|------------------|------------|-----|
| Conc. pg/mL | ECL Signal | %CV |
| 1890 | 311,000 | 4.7 |
| 473 | 82,400 | 4.3 |
| 118 | 20,700 | 5.8 |
| 29.5 | 5,260 | 5.2 |
| 7.38 | 1,460 | 5.5 |
| 1.85 | 411 | 6.1 |
| 0.461 | 147 | 4.0 |
| 0.00 | 58 | 7.5 |

CERTIFICATE OF ANALYSIS

Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|--------------------------------------------|------------------|---------------------|-----------------|
| Human TARC Plate | Z0046465 | 2-8°C | 30 Sep 2019 |
| Anti-hu TARC Antibody | D0080693 | 2-8°C | 30 Jun 2018 |
| Chemokine Panel 1 (human) Calibrator Blend | A0080122 | 2-8°C | 28 Feb 2022 |
| Diluent 43 | M0430017 | ≤ -10°C | 31 Dec 2018 |
| Diluent 3 | M0030070 | ≤ -10°C | 31 Jul 2018 |
| MSD Read Buffer T | Not Kit Specific | Not Kit Specific | N/A |

| Additional Components Included in V-PLEX Plus | |
|-----------------------------------------------|------------------|
| Plate Seals | Not Kit Specific |
| Wash Buffer (20x) | Not Kit Specific |
| Chemokine Panel 1 (human) Control 1 | Not Kit Specific |
| Chemokine Panel 1 (human) Control 2 | Not Kit Specific |
| Chemokine Panel 1 (human) Control 3 | Not Kit Specific |

Additional Comments:

| Assay | Spot | Specification | Comment |
|-----------|------|---------------|------------------------------------------------------------|
| 1 hu TARC | 1 | Control Conc. | Control 1 = 120.6%; Control 2 = 123.2%; Control 3 = 120.4% |

Statement:

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

| | Name | Function | Signature | Date |
|------------------------|---------------------|-------------------|----------------------------|-----------|
| Review/Approval | Vanessa Castleberry | Quality Assurance | <i>Vanessa Castleberry</i> | 13SEP2017 |