

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-Plex Plus Human IL-12p70 Kit
Kit Lot Number: K0081645
Expiration Date: 31 Oct 2022
Kit Catalog Numbers: K151QV-Series

Precision and Accuracy Specifications:

| Spot | Name | Calibrator | | Proinflammatory Panel 1 (human) Control 1 | | Proinflammatory Panel 1 (human) Control 2 | | Proinflammatory Panel 1 (human) Control 3 | |
|------|----------|------------|-------|---|-------|---|-------|---|-------|
| | | Recovery | CV | Recovery | CV | Recovery | CV | Recovery | CV |
| 1 | IL-12p70 | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% | 80% - 120% | < 20% |

| Spot | Name | LLOQ | | | ULOQ | | |
|------|----------|-------|------------|-------|-------|------------|-------|
| | | Value | Recovery | CV | Value | Recovery | CV |
| 1 | IL-12p70 | 1.22 | 80% - 120% | < 20% | 315 | 80% - 120% | < 20% |

Stock Concentration of Calibrator:


| Spot | Name | Conc. | Units |
|------|----------|-------|-------|
| 1 | IL-12p70 | 457 | pg/mL |

Lot-Specific Components:

| Description | Lot Number | Storage Temperature | Expiration Date |
|--|------------|---------------------|-----------------|
| Human IL-12p70 Plate | Z0047463 | 2-8°C | 31 Jul 2023 |
| Anti-hu IL-12p70 Antibody | D0081339 | 2-8°C | 31 Jan 2023 |
| Proinflammatory Panel 1 (human) Calibrator blend | A0080288 | 2-8°C | 31 Mar 2026 |
| Diluent 2 | M0020115 | ≤ -10°C | 31 May 2023 |
| Diluent 3 | M0030100 | ≤ -10°C | 31 Oct 2022 |

Additional components may be provided to execute the assay. Additional components are not lot-specific.

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

| | Name | Function | Signature | Date |
|-----------------|---------------|----------|--|-------------|
| Review/Approval | Karen Hamilla | Quality |  | 21 JUN 2021 |