

Certificate of Analysis

VeriPlex™ Human Interferon 9-Plex ELISA Kit

Catalog No: 51500-1

Lot No: 7302R

Expiration: December 31, 2021

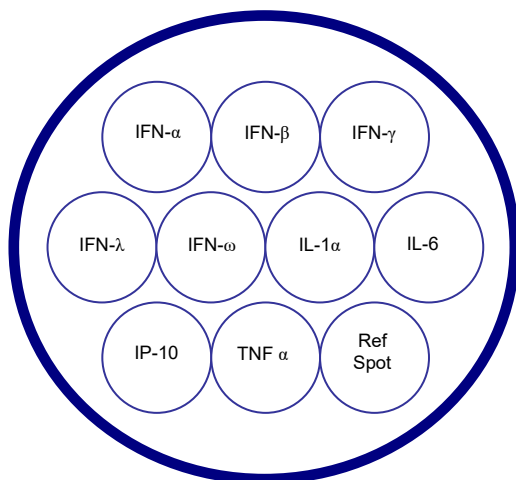
Software Code (input as product number in Q-View™ software): PCHIM180801-PB

| Kit Components | Part No. | Lot No. | Quantity | Reconstitution Instructions |
|--------------------------------------|----------|-------------|----------|---|
| Plate(s) | 311952HU | PCHIP201023 | 1 | N/A |
| Plate Sealers | N/A | N/A | 3 | N/A |
| Wash Solution Concentrate | 101158GR | PCHW200813 | 1 bottle | 950 ml DI Water |
| Human IFN Multiplex Antigen Standard | 311955HU | PCHIM180801 | 1 vial | 0.600 ml Sample Diluent |
| Human IFN Multiplex Detection Mix | 311956HU | PCHID201002 | 1 vial | 6 ml DI Water |
| HRP Concentrate | 101173GR | PCHN200707 | 6 ml | N/A |
| Sample Diluent | 311969HU | PCHIS200319 | 10 ml | N/A |
| Substrate A | 101159GR | PCHA201007 | 3 ml | Mix equal parts of A and B together and store in the dark |
| Substrate B | 101160GR | PCHB201007 | 3 ml | |
| Diluent Additive II | SMP208-1 | DA075 | 1 vial | N/A |
| Diluent Additive III | SMP209-3 | DA076 | 1 vial | N/A |

| Kit Components | Part No. | Lot No. | Quantity | Reconstitution Instructions |
|----------------|----------|------------|----------|--|
| Assay Diluent | 311983HU | PCHV200702 | 1 vial | 5.0 ml DI Water + 0.480 ml Diluent Additive II + 0.471 ml Diluent Additive III to reconstitute Assay Diluent |

Spot Location in each well:

Assay Range for Analytes:



| Analyte | Assigned High Value (pg/ml) | Assay Range (pg/ml) | LLOD (pg/ml) | LLOQ (pg/ml) |
|---------|-----------------------------|---------------------|--------------|--------------|
| IFN-α | 3400 | 3400 – 4.7 | 0.8 | 12.2 |
| IFN-β | 20100 | 20100 – 27.6 | 5.6 | 60.1 |
| IFN-γ | 5600 | 5600 – 7.7 | 3.2 | 8.9 |
| IFN-λ | 12900 | 12900 – 17.7 | 3.4 | 35.2 |
| IFN-ω | 3800 | 3800 – 5.2 | 0.9 | 5.2 |
| IL-1α | 1100 | 1100 – 1.5 | 0.3 | 1.5 |
| IL-6 | 22000 | 22000 – 30.2 | 3.3 | 30.2 |
| IP-10 | 1400 | 1400 – 1.9 | 0.7 | 21.1 |
| TNF-α | 2900 | 2900 – 4.0 | 0.0 | 3.7 |

Authorization

Released by: _____

Date: May 10, 2021

Sold under license from Pestka Biomedical Laboratories, Inc. d/b/a PBL Assay Science. For research use only. Not for diagnostic or clinical use in, or administration to, humans. Not for resale in original or any modified form, including inclusion in a kit, for any purpose. Not for use in the preparation of any commercial product.