



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA- 07418

Product Description: U-PLEX[®] Human IL-22 Antibody Set

Catalog Numbers: B21WI-2, B21WI-3

Antibody Set Lot Number: SET00134A

Expiration Date: 28 FEB 2019

| U-PLEX [®] Antibody Set Components | | | |
|---|------------|---------------------|-----------------|
| Description | Lot Number | Storage Temperature | Expiration Date |
| Biotin Anti-hu IL-22 Antibody | B00U0134A | 2-8°C | 28 FEB 2019 |
| SULFO-TAG [™] Anti-hu IL-22 Antibody | D00U0134A | 2-8°C | 28 FEB 2019 |

Statement: The above antibody set was tested using MSD procedures and conforms to MSD's specifications. It 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> | 14 OCT 2016 |