



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-07275

Product Description: U-PLEX[®] Mouse IL-21 Antibody Set

Catalog Numbers: B22WB-2, B22WB-3

Antibody Set Lot Number: SET00130A

Expiration Date: 31 JAN 2019

U-PLEX [®] Antibody Set Components			
Description	Lot Number	Storage Temperature	Expiration Date
Biotin Anti-ms IL-21 Antibody	B00U0130A	2-8°C	31 JAN 2019
SULFO-TAG [™] Anti-ms IL-21 Antibody	D00U0130A	2-8°C	31 JAN 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>	06 SEP 2016