



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-05631

Product Description: U-PLEX[®] Mouse IL-6 Antibody Set


Catalog Numbers: B22TX-2, B22TX-3

Antibody Set Lot Number: SET00052A

Expiration Date: 31 OCT 2017

U-PLEX [®] Antibody Set Components			
Description	Lot Number	Storage Temperature	Expiration Date
Biotin Anti-ms IL-6 Antibody	B00U0052A	2-8°C	31 OCT 2017
SULFO-TAG [™] Anti-ms IL-6 Antibody	D00U0052A	2-8°C	31 OCT 2017

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	Anastasia Belilovets	Quality Assurance		21 Jul 2015