



## CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-06732

**Product Description:** U-PLEX<sup>®</sup> Mouse IL-17E/IL-25 Antibody Set

**Catalog Numbers:** B22VZ-2, B22VZ-3

**Antibody Set Lot Number:** SET00106A

**Expiration Date:** 31 JUL 2018

U-PLEX <sup>®</sup> Antibody Set Components			
Description	Lot Number	Storage Temperature	Expiration Date
Biotin Anti-ms IL-17E/IL-25 Antibody	B00U0106A	2-8°C	31 AUG 2018
SULFO-TAG <sup>®</sup> Anti-ms IL-17E/IL-25 Antibody	D00U0106A	2-8°C	31 JUL 2018

**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>	02 MAY 2016