



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-06562

Product Description: U-PLEX[®] Human IL-17E/IL-25 Antibody Set

Catalog Numbers: B21VZ-2, B21VZ-3

Antibody Set Lot Number: SET00090A

Expiration Date: 31 JUL 2018

U-PLEX [®] Antibody Set Components			
Description	Lot Number	Storage Temperature	Expiration Date
Biotin Anti-hu IL-17E/IL-25 Antibody	B00U0090A	2-8°C	31 JUL 2018
SULFO-TAG [®] Anti-hu IL-17E/IL-25 Antibody	D00U0090A	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>	06 MAY 2016