



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA- 07415

Product Description: U-PLEX[®] Human MIP-3 β Antibody Set

Catalog Numbers: B21VA-2, B21VA-3

Antibody Set Lot Number: SET00133A

Expiration Date: 28 FEB 2019

U-PLEX [®] Antibody Set Components			
Description	Lot Number	Storage Temperature	Expiration Date
Biotin Anti-hu MIP-3 β Antibody	B00U0133A	2-8°C	28 FEB 2019
SULFO-TAG [™] Anti-hu MIP-3 β Antibody	D00U0133A	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