



## CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-07222

**Product Description:** U-PLEX<sup>®</sup> Mouse GM-CSF Antibody Set

**Catalog Numbers:** B22UM-2, B22UM-3

**Antibody Set Lot Number:** SET00132A

**Expiration Date:** 31 JAN 2019

U-PLEX <sup>®</sup> Antibody Set Components			
Description	Lot Number	Storage Temperature	Expiration Date
Biotin Anti-ms GM-CSF Antibody	B00U0132A	2-8°C	31 JAN 2019
SULFO-TAG <sup>™</sup> Anti-ms GM-CSF Antibody	D00U0132A	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>	17AUG2016