



CERTIFICATE OF ANALYSIS

COA - 07656

Product Description: SULFO-TAG™ Anti-Hu/NHP IgG Antibody

Catalog #/Size: D20JL-6/100 µg

Lot Number: W0016710S

Expiration Date: 31 MAR 2019

Statement: This product was manufactured and tested according to approved MSD documents. This lot meets MSD's specifications.

Volume: 200 µL



Concentration: 500 µg/mL

L/P: 3.0

Diluent: Diluent 100 w/ Azide

Storage conditions: 2-8°C

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Preparer	Meredith Hutchinson	Production Control		13 JAN 2017
Review/Approval	Dawn Stollar	Quality Assurance		17 JAN 2017