



## CERTIFICATE OF ANALYSIS

COA - 05263

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

**Catalog #/Size:** D20JL-6/100 µg

**Lot Number:** W0015481S

**Expiration Date:** 30 APR 2017

**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

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 Yeager	Inventory Technician		06 MAR 2015
Review/Approval	Staci Leffel	Manufacturing Manager		10 MAR 2015