



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: SULFO-TAG™ Anti-Hu/NHP Kappa Light Chain Antibody

Catalog #/Size: D20TF -Series

Lot Number: W0017251S

Expiration Date: 30 NOV 2019

Statement: This product was manufactured and tested according to approved MSD documents.

Total Volume: 200 µL

Concentration: 500 µg/mL

L/P: 8.7

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
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	09 OCT 2017