



CERTIFICATE OF ANALYSIS

COA - 04742

Product Description: MSD SULFO-TAG™ NHS Ester

Catalog #/Size: R91AN-1 / 150 nmol

Lot Number: S0080058



Expiration Date: 31 JUL 2016

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

Purity (by HPLC): > 90%

Quantity: 100% to 120% of nominal

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

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biologics/Chemistry		11 SEP 2014
Review/Approval	Anastasia Belilovets	Quality Control		11 SEP 2014