



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD® TAG NHS-Ester

Catalog #/Size: R91BN-2 / 500 nmol

Lot Number: T0040009

Expiration Date: 31 AUG 2026

Storage Temperature: $\leq -10^{\circ}\text{C}$

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


Purity specification (by HPLC): $> 90\%$

Measured Average Purity: 99%

Quantity: 100% to 120% of nominal

Measured Average Quantity: 113%

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		03 AUG 2021