



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-07286

Product Description: U-PLEX[®] 10-Assay, 96-Well SECTOR[®] Plate
Catalog Number: N05235A-1
Lot Number: Z00U0023
Expiration Date: 31 JAN 2019
Storage Condition: 2-8°C

Acceptance Data:

Metric	Specifications	All Spots
Average Intraplate CV	≤ 10.0%	Pass
Average Interplate CV	≤ 13.0%	Pass
Signal Range at Binding Capacity	≤ 30.0%	Pass

Statement: The above plate lot was tested using MSD procedures and conforms to MSD's specifications. It 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>	12 SEP 2016