

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: U-PLEX® 8-Assay, 96-Well SECTOR® Plate **Catalog Number:** N05233A-1 **Lot Number:** Z00U0051G **Expiration Date:** 30 APR 2020 **Storage Condition:** 2-8°C

Acceptance Data:

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

Statement: The above plate lot is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	Dawn Stollar	01 DEC 2017