

**CERTIFICATE OF ANALYSIS**

Product Description: T1D Autoantibody Panel 1 (hu) 96-Well 10-Spot SECTOR Plate
Catalog Number: N05757A-1
Lot Number: Z0048195
Expiration Date: 30 APR 2028
Storage Temperature: 2-8°C

Acceptance Data:

Metric	Specifications	All Spots
Average Intraplate CV	$\leq 10.0\%$	Pass
CV of Intraplate Averages	$\leq 13.0\%$	Pass

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		13 JAN 2026