

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX T1D Autoantibody Panel 1 (human) Kit, QuickPlex**Lot Number:** K0082877**Expiration Date:** 31 MAR 2028**Catalog Number:** K15757-Series**Lot-Specific Components:**

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
T1D Autoantibody Panel 1 (hu) 96-Well 10-Spot QuickPlex Ultra Plate	Z0048196	2-8°C	30 APR 2028
Biotin Human GAD2/GAD-65	565369	≤ -70°C	30 JUN 2028
Biotin Human ZnT8	565376	≤ -70°C	30 JUN 2028
Biotin Human IA-2/R-PTP-N	565380	≤ -70°C	30 JUN 2028
Biotin Human Insulin	565383	≤ -70°C	30 JUN 2028
SULFO-TAG Human GAD2/GAD-65	D0082561	≤ -70°C	31 MAY 2028
SULFO-TAG Human ZnT8	D0082563	≤ -70°C	31 MAY 2028
SULFO-TAG Human IA-2/R-PTP-N	D0082565	≤ -70°C	31 MAY 2028
SULFO-TAG Human Insulin	D0082567	≤ -70°C	31 MAY 2028
T1D Autoantibody Panel 1 (human) Calibrator Blend	A0080447	≤ -70°C	30 APR 2028
Diluent 6	M0060047	≤ -10°C	31 MAR 2028

Additional components may be provided to execute the assay. Additional components are not lot-specific.

Precision, Accuracy, and Sensitivity Specifications:

Spot	Name	Calibrator		T1D Autoantibody Panel 1 (human) Control 1		T1D Autoantibody Panel 1 (human) Control 2	
		Recovery	CV	Recovery	CV	Recovery	CV
1	GADA	80% - 120%	< 20%	70% - 130%	< 20%	70% - 130%	< 20%
3	IA-2A	80% - 120%	< 20%	70% - 130%	< 20%	70% - 130%	< 20%
8	ZnT8A	80% - 120%	< 20%	70% - 130%	< 20%	70% - 130%	< 20%
10	IAA	80% - 120%	< 20%	70% - 130%	< 20%	70% - 130%	< 20%



CERTIFICATE OF ANALYSIS

Spot	Name	LLOQ			ULOQ			LLOD
		Value	Recovery	CV	Value	Recovery	CV	Value
1	GADA	9.00	75% - 125%	<20%	2000	75% - 125%	<20%	3.00
3	IA-2A	7.00	75% - 125%	<20%	900	75% - 125%	<20%	4.00
8	ZnT8A	16.0	75% - 125%	<20%	1000	75% - 125%	<20%	8.00
10	IAA	18.0	75% - 125%	<20%	1200	75% - 125%	<20%	12.0

Non-Specific Binding Specifications:

Spot	Name	Spec
1	GADA	< 1.0%
3	IA-2A	< 1.0%
8	ZnT8A	< 1.0%
10	IAA	< 1.0%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	GADA	25200	U/mL
3	IA-2A	11400	U/mL
8	ZnT8A	14200	U/mL
10	IAA	23450	U/mL

Dilution to top of Curve: 5X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		25 MAR 2026