

**CERTIFICATE OF ANALYSIS**

Description: T1D Autoantibody Panel 1 (human) Controls
Catalog Number: C5757-1, C5757-2
Lot Number: A00C1034, A00C1035
Expiration Date: 31 MAY 2028

Components:

Description	Lot Number	Storage Temperature	Expiration Date
T1D Autoantibody Panel 1 (human) Control 1	A00C1034	$\leq -70^{\circ}\text{C}$	31 MAY 2028
T1D Autoantibody Panel 1 (human) Control 2	A00C1035	$\leq -70^{\circ}\text{C}$	31 MAY 2028

Concentration Assignment:

Analyte	T1D Autoantibody Panel 1 (human) Control 1	T1D Autoantibody Panel 1 (human) Control 2	Unit of Measure
GADA	707	137	U/mL
IA-2A	320	52.8	U/mL
ZnT8A	629	158	U/mL
IAA	533	172	U/mL

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		27 JAN 2026