



## CERTIFICATE OF ANALYSIS

Page 1 of 0

**Product Description:** Human Type 1 Diabetes Antibody Panel 1 Control Pack

**Catalog Number:** C01XX-2

**Lot Number:** CP000095

**Expiration Date:** 31 MAY 2022

**Safety Consideration:** Biosafety level 2 (BSL-2) material; handle according to appropriate safety considerations.

### Concentration of Materials Tested:

Units: U/mL	Assay				
	Anti-TGM2	Anti-GAD65	Anti-ZnT8	Anti-Insulin	Anti-IA-2
Positive Control 1 (A00C0694)	86.4	901	114	23.4	9.77
Positive Control 2 (A00C0695)	33.7	143	36.9	3.32	4.44
Negative Control (A00C0696)	< 12.0	< 14.0	< 6.00	< 1.50	< 1.50

Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Human T1D Antibody Panel 1 Positive Control 1	A00C0694	≤ -70°C	31 MAY 2022
Human T1D Antibody Panel 1 Positive Control 2	A00C0695	≤ -70°C	31 MAY 2022
Human T1D Antibody Panel 1 Negative Control	A00C0696	≤ -70°C	31 MAY 2022

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		20 NOV 2019