

**\*Confidential – Akston Biosciences**  
**Cross-Reactivity Panel Results – Final**  
**Run at Tier 1 University Hospital System**

AntiCoV-ID™ IgG ELISA - Akston			
Serum Panel Containing IgG's Against	N	Negative	%Negative
Adenovirus E	4	4	100%
Coronavirus 229E	8	8	100%
Coronavirus HKU1	6	6	100%
Coronavirus NL63	9	8	89%
Coronavirus OC43	15	15	100%
Coxsackie/Echovirus	12	12	100%
Influenza A	2	2	100%
Influenza B	6	6	100%
Influenza A, novel H1N1	8	8	100%
Parainfluenzavirus 1	2	2	100%
Parainfluenzavirus 3	2	2	100%
Parainfluenzavirus 4	2	2	100%
Metapneumovirus (A,B)	2	2	100%
Rhinovirus	11	11	100%
Rhinovirus & Coxsackie/Echovirus	14	14	100%
RSVA	2	2	100%
<b>Total</b>	<b>105</b>	<b>104</b>	
<b>False Positives</b>	<b>1</b>		
<b>False Positive Rate in Panel</b>	<b>&lt; 1.0%</b>		

EuroImmune IgG ELISA			
Serum Panel Containing IgG's Against	N	Negative	%Negative
Adenovirus E	4	4	100%
Coronavirus 229E	8	8	100%
Coronavirus HKU1	6	6	100%
Coronavirus NL63	9	9	100%
Coronavirus OC43	15	15	100%
Coxsackie/Echovirus	12	11	92%
Influenza A	2	2	100%
Influenza B	6	4	67%
Influenza A, novel H1N1	8	7	88%
Parainfluenzavirus 1	2	2	100%
Parainfluenzavirus 3	2	2	100%
Parainfluenzavirus 4	2	2	100%
Metapneumovirus (A,B)	2	2	100%
Rhinovirus	11	11	100%
Rhinovirus & Coxsackie/Echovirus	14	12	86%
RSVA	2	2	100%
<b>Total</b>	<b>105</b>	<b>99</b>	
<b>False Positives</b>	<b>6</b>		
<b>False Positive Rate in Panel</b>	<b>5.7%</b>		