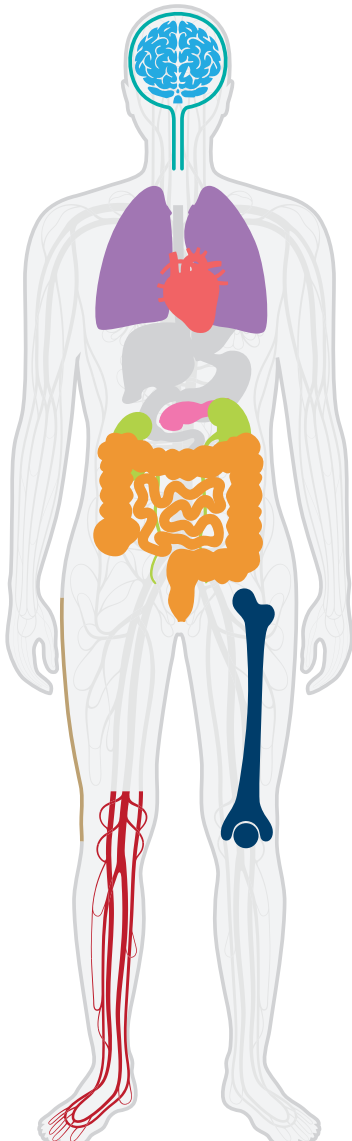




LONG COVID IMPACTS






Mental Health

- Anxiety
- Depression
- Sleep problems
- Substance abuse




Respiratory System

- Cough
- Low blood oxygen
- Shortness of breath




Kidney

- Acute kidney injury
- Chronic kidney disease




Gastrointestinal

- Diarrhea
- Acid reflux
- Constipation




Skin Disorders

- Rash
- Hair loss




Blood Disorders

- Anemia
- Blood clots




Nervous System

- Stroke
- Headaches
- Memory problems
- Loss of smell and taste




Cardiovascular

- Arrhythmia
- Palpitations
- Heart failure
- Acute coronary disease




Metabolic/Endocrine

- Obesity
- Diabetes
- High cholesterol



Musculoskeletal

- Joint pain
- Muscle weakness



General

- Fatigue
- Malaise
- Mitochondrial dysfunction

Use the links below to browse products and resources from Cayman related to Long COVID research.

ANXIETY	THROMBOSIS	OBESITY
DEPRESSION	ERYTHROCYTES & HEMOGLOBIN	DIABETES
SLEEP & BIOLOGICAL RHYTHMS	STROKE	DYSLIPIDEMIAS
ADDICTION RESEARCH	NEUROPROTECTION	CHOLESTEROL METABOLISM
PULMONARY DISEASES	LEARNING & MEMORY	PAIN RESEARCH
KIDNEY & RENAL DISEASE	CORONARY ARTERY DISEASE	MITOCHONDRIAL BIOLOGY
GASTRIC DISEASE	HEART FAILURE	CORONAVIRUS RESEARCH
COAGULATION & HEMOSTASIS	ARRHYTHMIA	