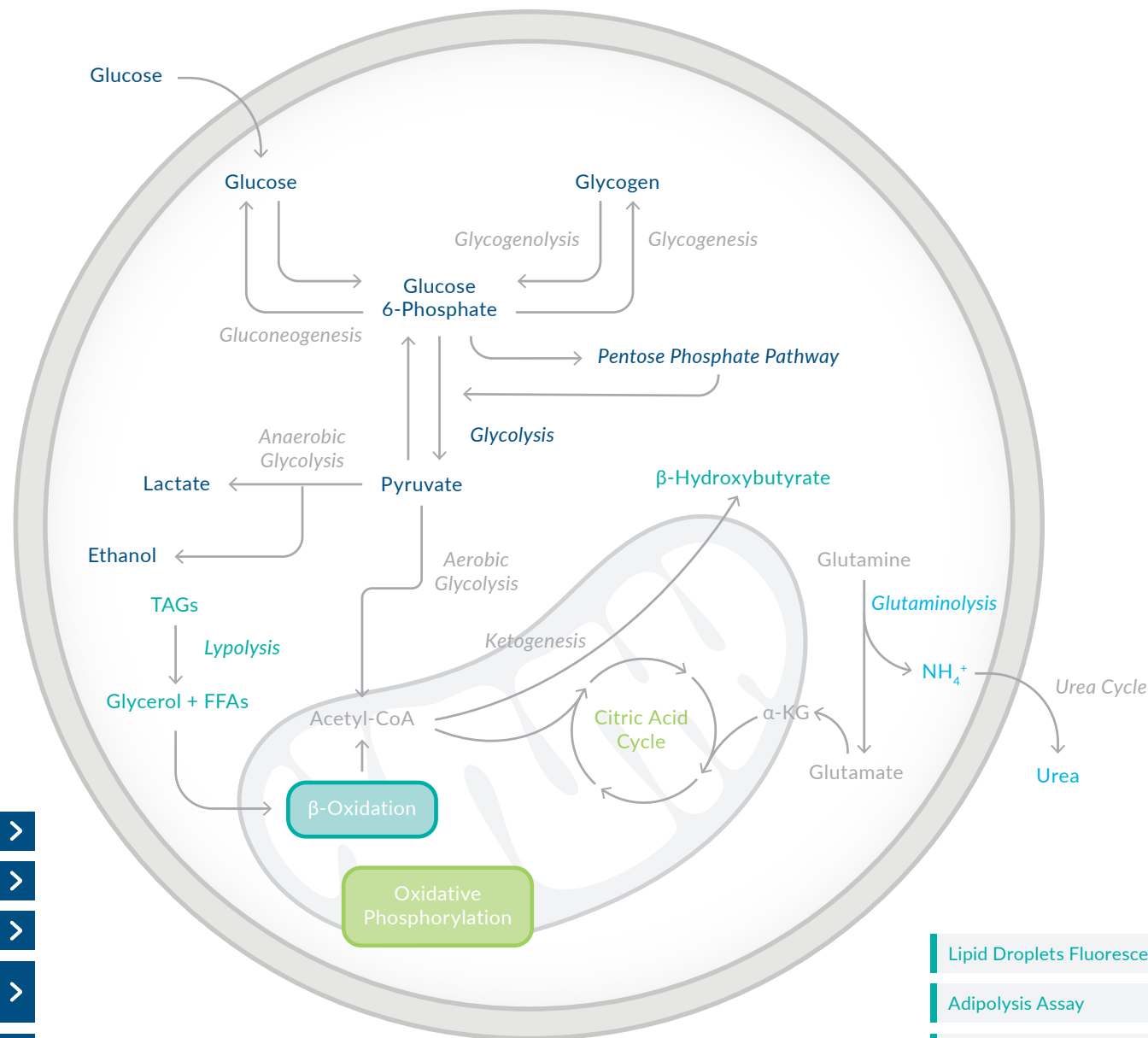


Cellular Metabolism Assay Kits

© 2025 Cayman Chemical



Category Key

= Carbohydrate Metabolism

= Amino Acid Metabolism

= Lipid Metabolism

= Aerobic Cellular Respiration

Glucose Uptake Cell-Based Assay >

Glucose Colorimetric Assay >

Glucose-6-Phosphate Fluorometric Assay >

Glucose-6-Phosphate Dehydrogenase Activity Assay >

Glycolysis Cell-Based Assay >

NAD/NADH Cell-Based Assay >

Pyruvate Assay >

Lactate Assays >

Ethanol Colorimetric Assay >

Glycogen Assay >

Glutaminolysis Assays >

Ammonia Colorimetric Assay >

Urea Fluorometric Assay >

ATP Detection Assay - Luminescence >

Extracellular Oxygen Consumption Assay >

MitoCheck® Complex Activity Assays >

α-Ketoglutarate Detection Assay >

Mitochondrial Membrane Potential Assays >

MitoCheck® Citrate Synthase Activity Assay >

Aconitase Assay >

Lipid Droplets Fluorescence Assay >

Adipolysis Assay >

Triglyceride Colorimetric Assay >

Lipase Activity Assay >

Glycerol Colorimetric Assay >

Glycerol Cell-Based Assay >

Free Fatty Acid Fluorometric Assay >

β-Hydroxybutyrate Assays >