

# Cellular Energy Metabolism



G6P = glucose-6-phosphate  
 αKG = α-ketoglutarate  
 Ac-CoA = acetyl coenzyme A

A better way to *explore*

Learn more at [promega.com/EnergyMetabolism](http://promega.com/EnergyMetabolism)