

Combining Seahorse XF analysis with stable isotope tracing to reveal novel drug targets for metabolic and neurodegenerative disease

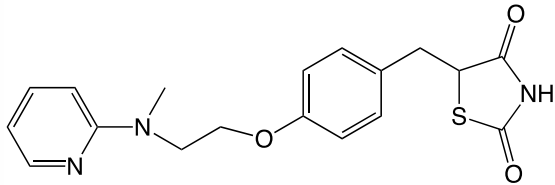
Ajit Divakaruni, Ph.D.

July 20, 2017

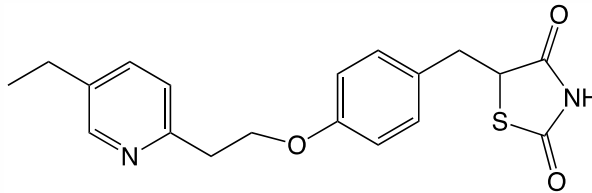
Asst. Professor of Molecular & Medical Pharmacology, UCLA

Thiazolidinediones (TZDs)

Rosiglitazone (Avandia®)



Pioglitazone (Actos®)

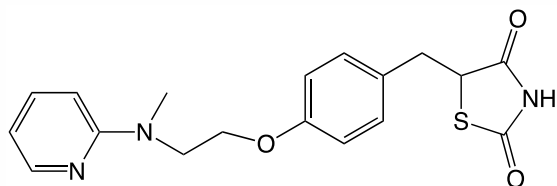


Effective insulin sensitizers with prohibitive side effects

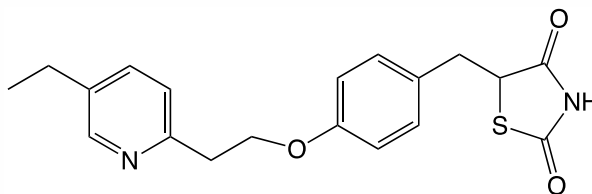
Is PPAR γ agonism the only mechanism of action?

Thiazolidinediones (TZDs)

Rosiglitazone (Avandia®)



Pioglitazone (Actos®)

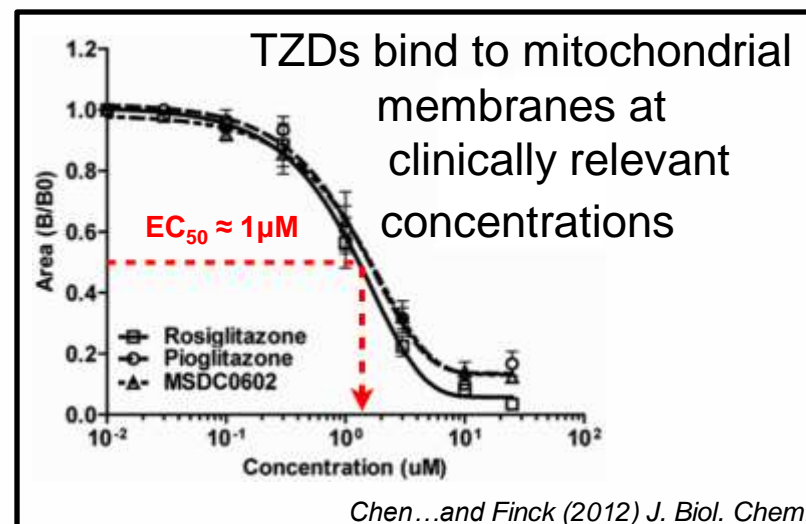


Effective insulin sensitizers with prohibitive side effects

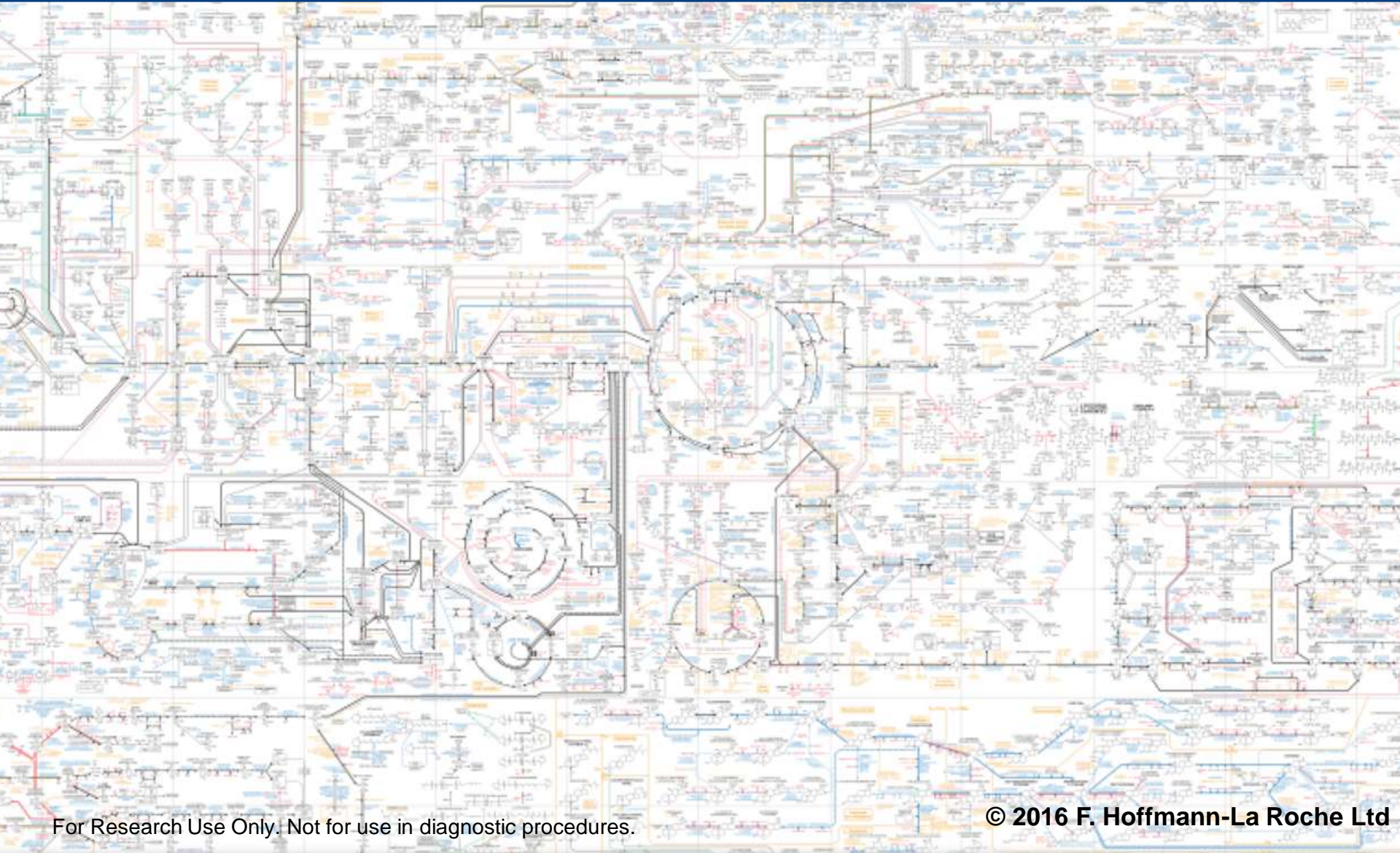
Is PPAR γ agonism the only mechanism of action?

Rapid *in vitro* and *in vivo* effects

An insulin-sensitizing effect persists
in some tissue-specific KO models



Metabolism



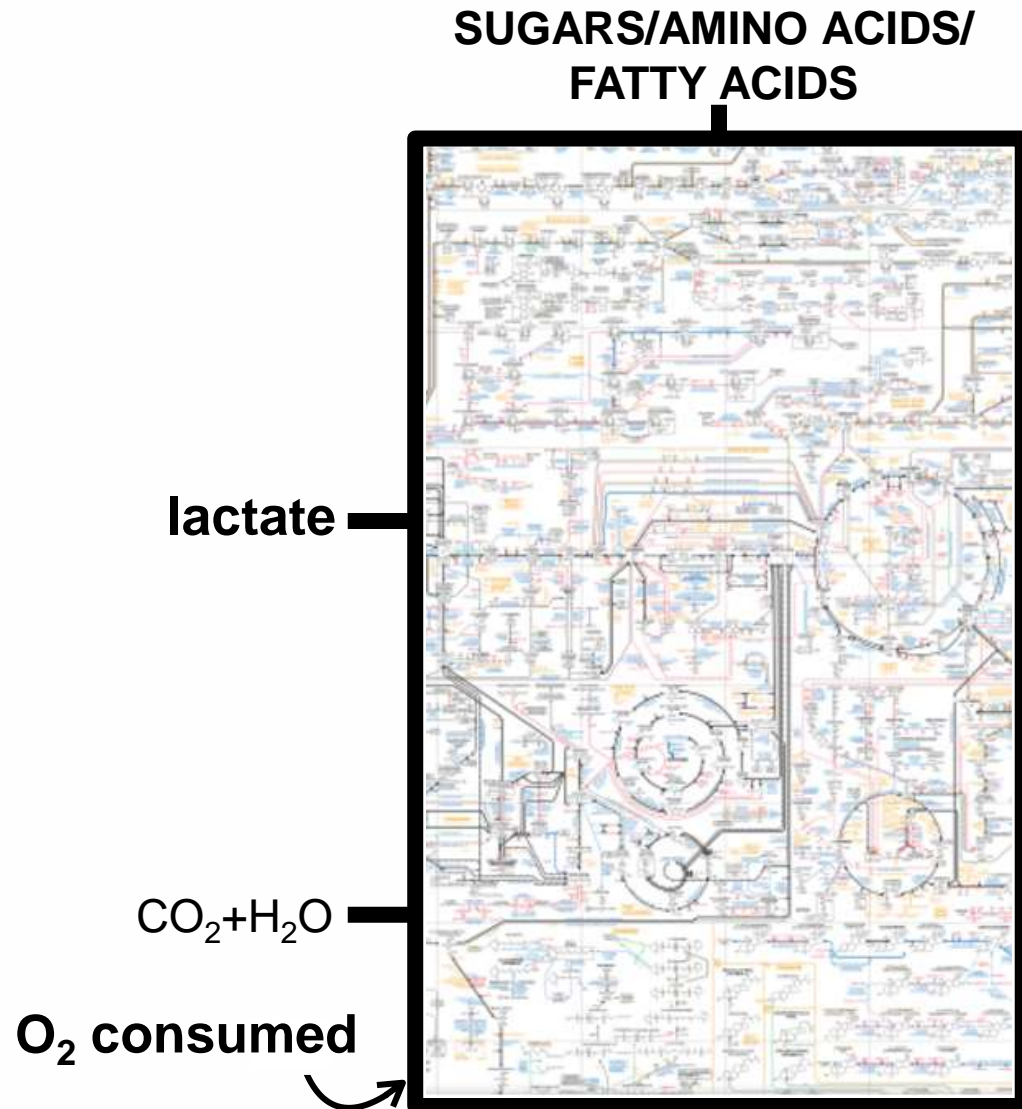
A “black box” approach

(1) MEASURE RATES OF INPUTS AND OUTPUTS

(Respiration, lactate efflux)

Quantitative flux through
an intact network

Real-time rates upon
acute perturbations



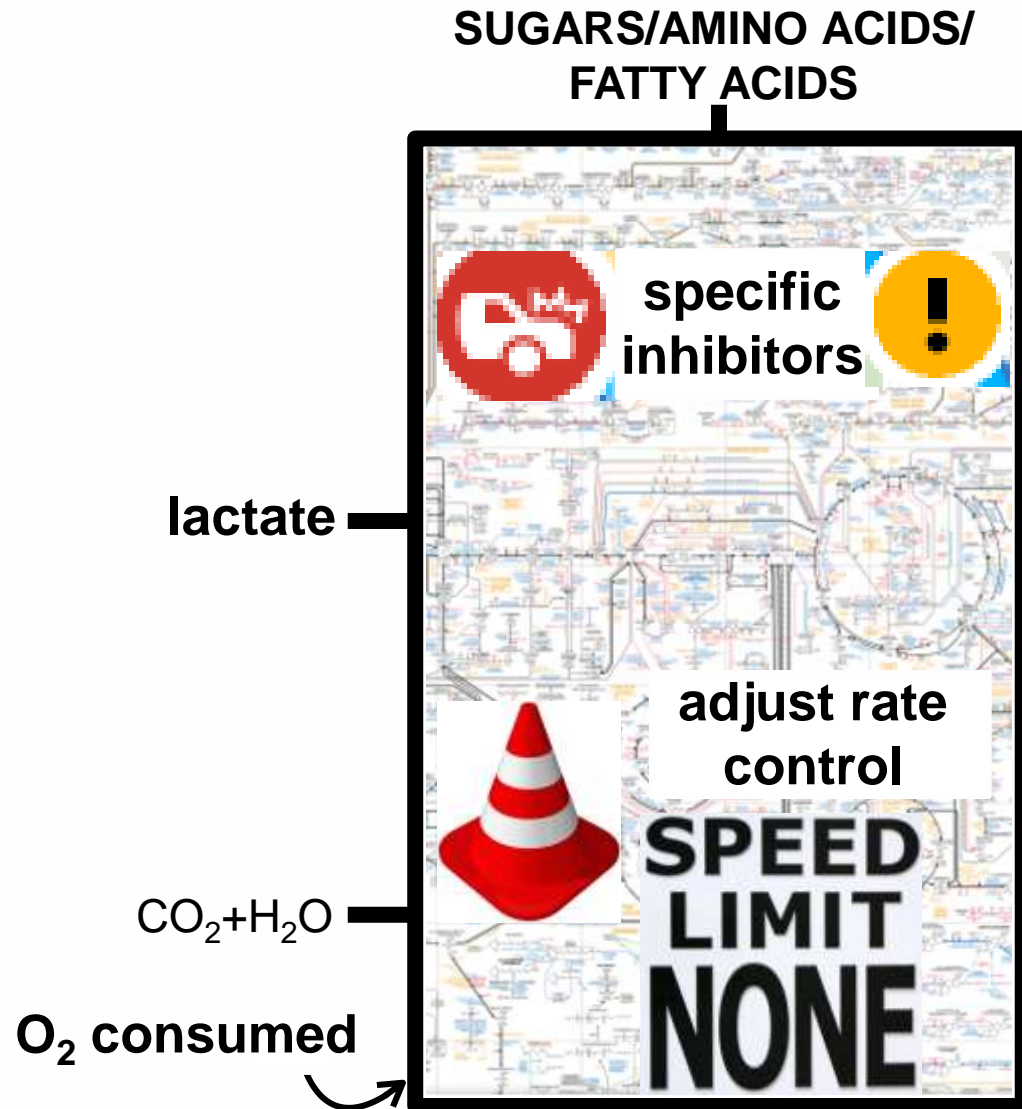
A “black box” approach

(1) MEASURE RATES OF INPUTS AND OUTPUTS

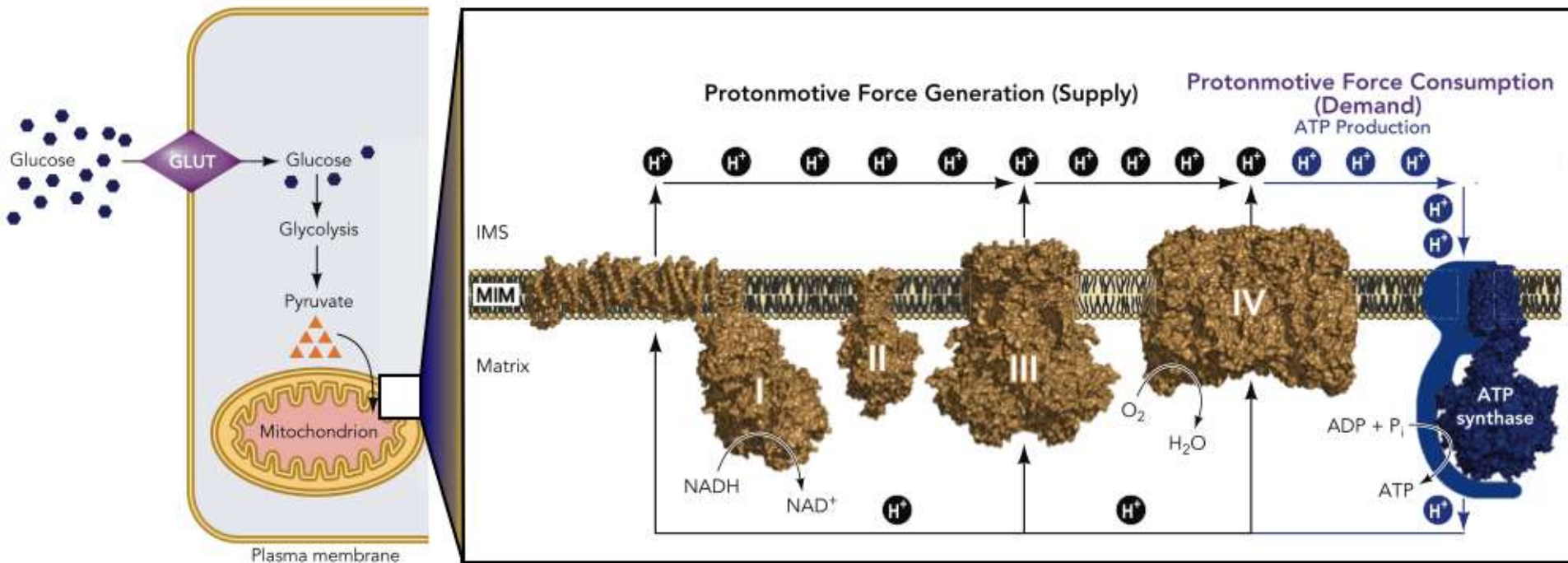
(Respiration, lactate efflux)

Quantitative flux through
an intact network

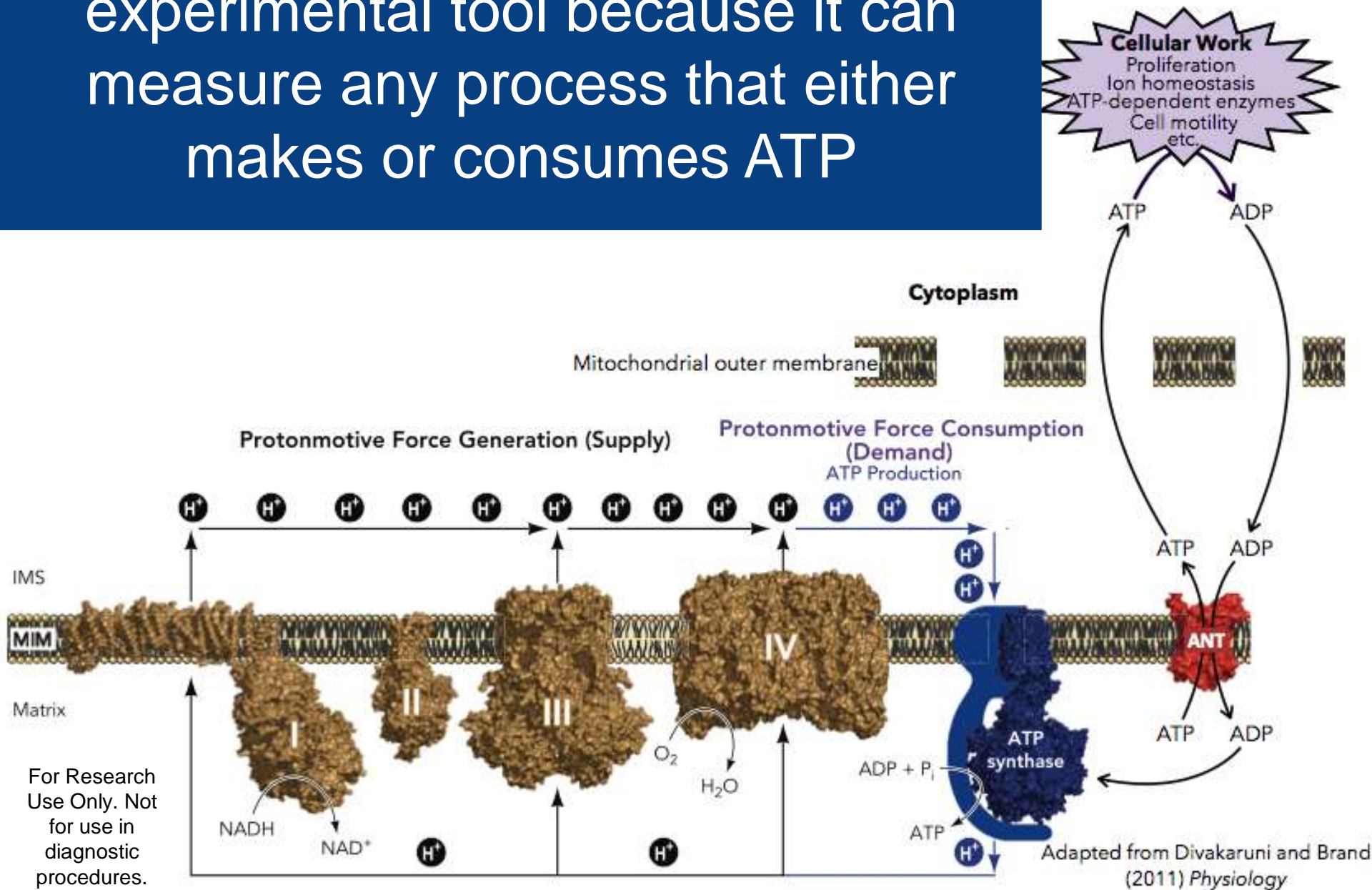
Real-time rates upon
acute perturbations



Respiration is a powerful experimental tool to study mitochondrial metabolism

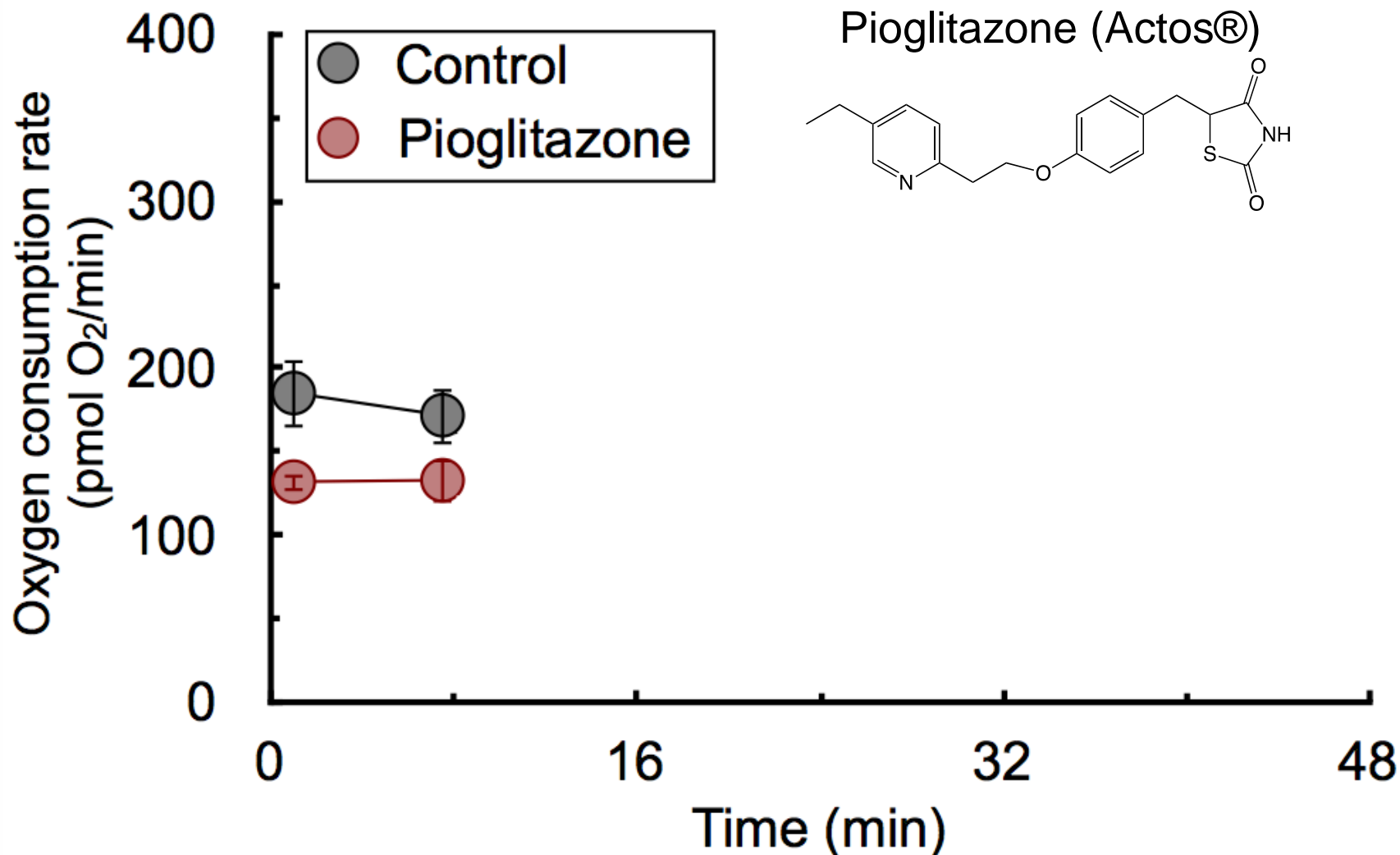


Respiration is a powerful experimental tool because it can measure any process that either makes or consumes ATP



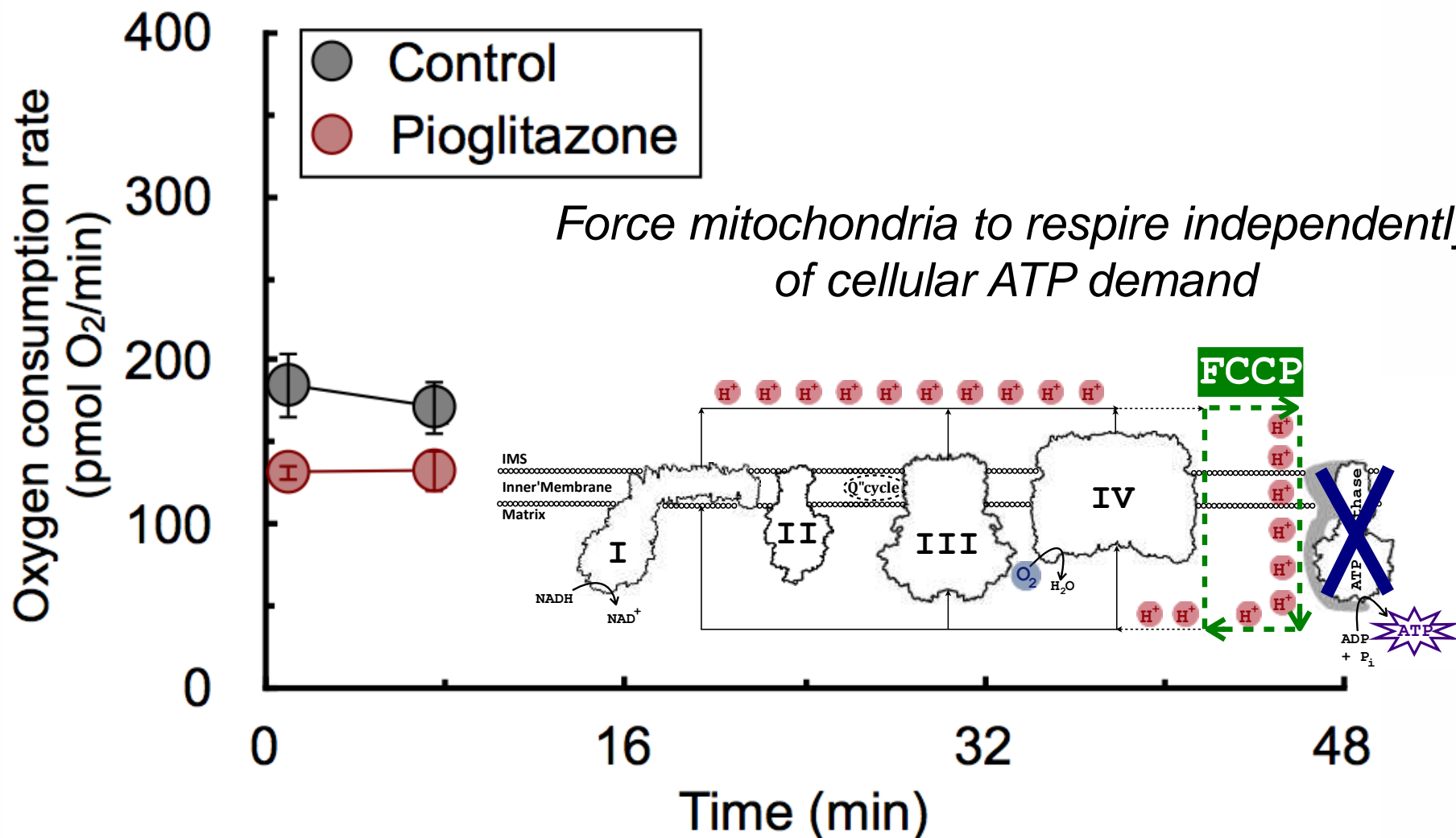
Is there a direct effect of TZDs on mitochondrial function?

RESPIRATION IN CULTURED MYOCYTES

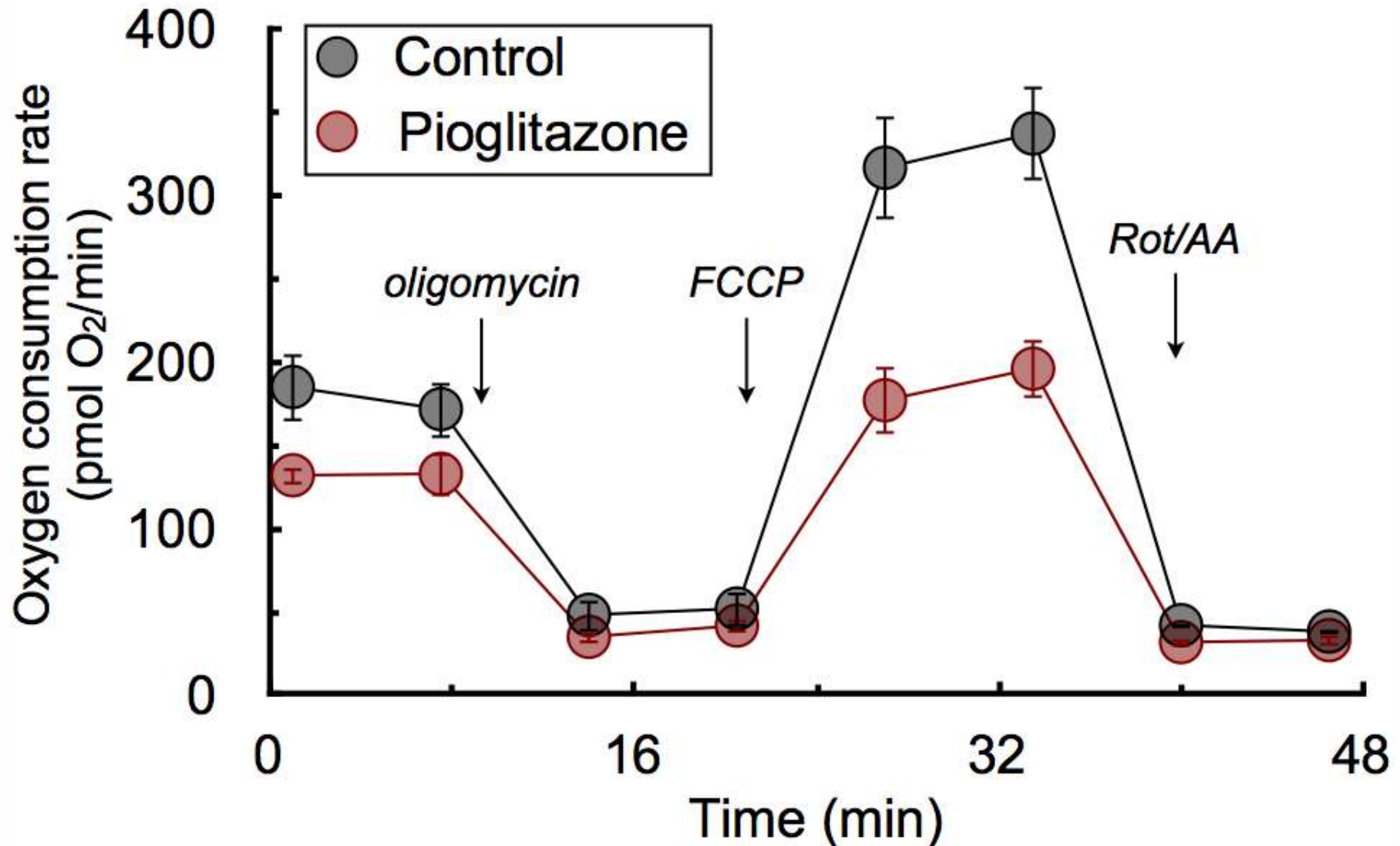


Is there a direct effect of TZDs on mitochondrial function?

RESPIRATION IN CULTURED MYOCYTES



Pioglitazone has a direct, acute effect on mitochondrial function



Decreased respiratory capacity could be due to dozens of possibilities

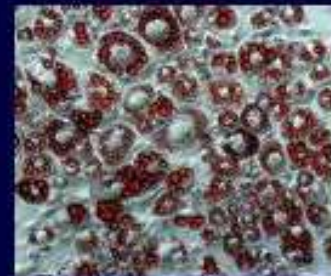
•SUBSTRATE/ION TRANSPORTERS

- GLUT •ARALAR •MCU •MCT •Plasma mem./mito. amino acid transport •CACT
- Mito. pyruvate carrier •Mito. dicarboxylate carrier •NCX •Mal./asp. shuttle •CD36



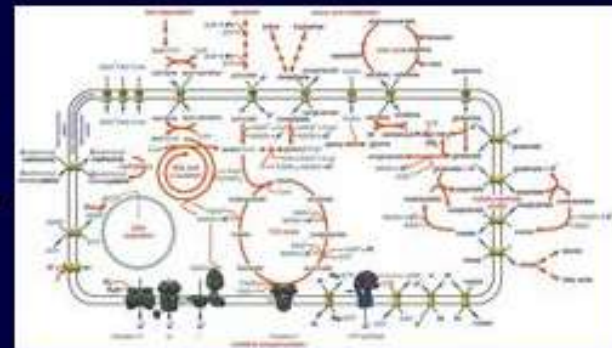
•MOBILIZATION OF INTERNAL ENERGY STORES

- Glycogenolysis •Endogenous lipolysis •Oxidation of endogenous amino acids



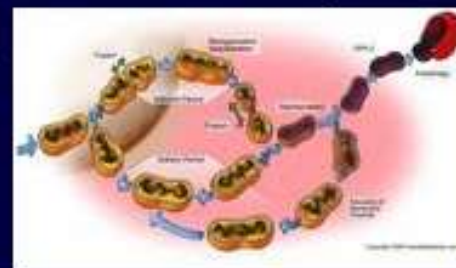
•ACTIVITY OF RATE-CONTROLLING ENZYME

- Respiratory complex I •PDH •CPT-1 •Hexokinase • α KGDH • β -oxidation enzymes
- ETFDH •Respiratory complex III •BCKDH •SDH •Aconitase •MDH •FH •CPT-2
- Ketone body oxidation enzymes •GAPDH •Citrate synthase •Cytochrome oxidase
- IDH •Glutaminase •GDH •Cytochrome c •AsAT •G3PDH •AIAT •BCAT



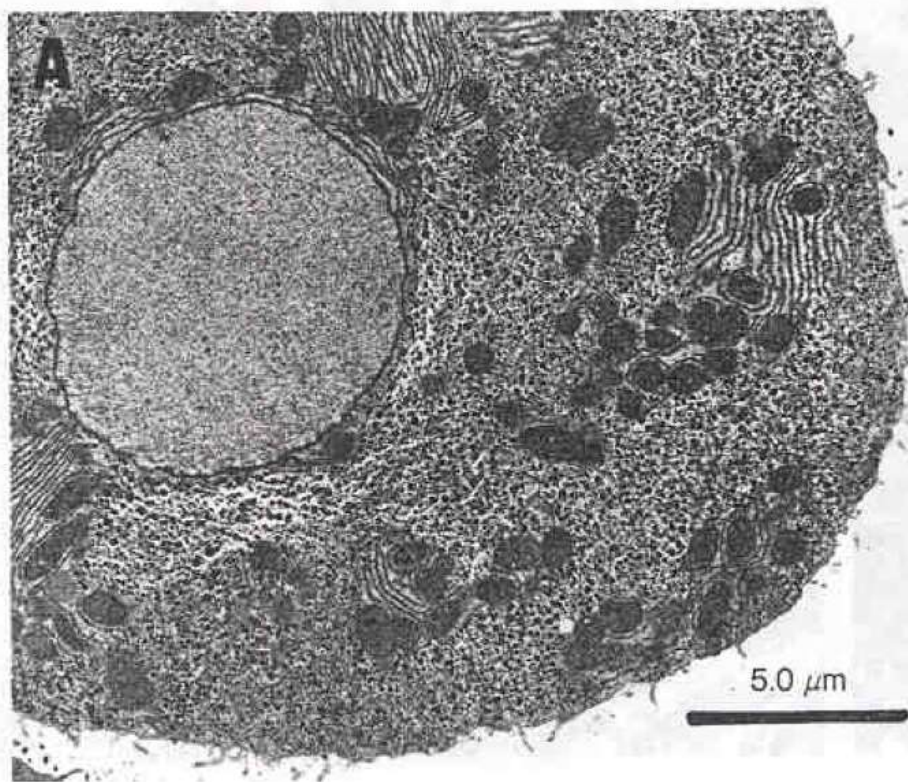
•MITOCHONDRIAL ULTRASTRUCTURE

- Content •Volume •Cristae density •Turnover •Fusion/Fission

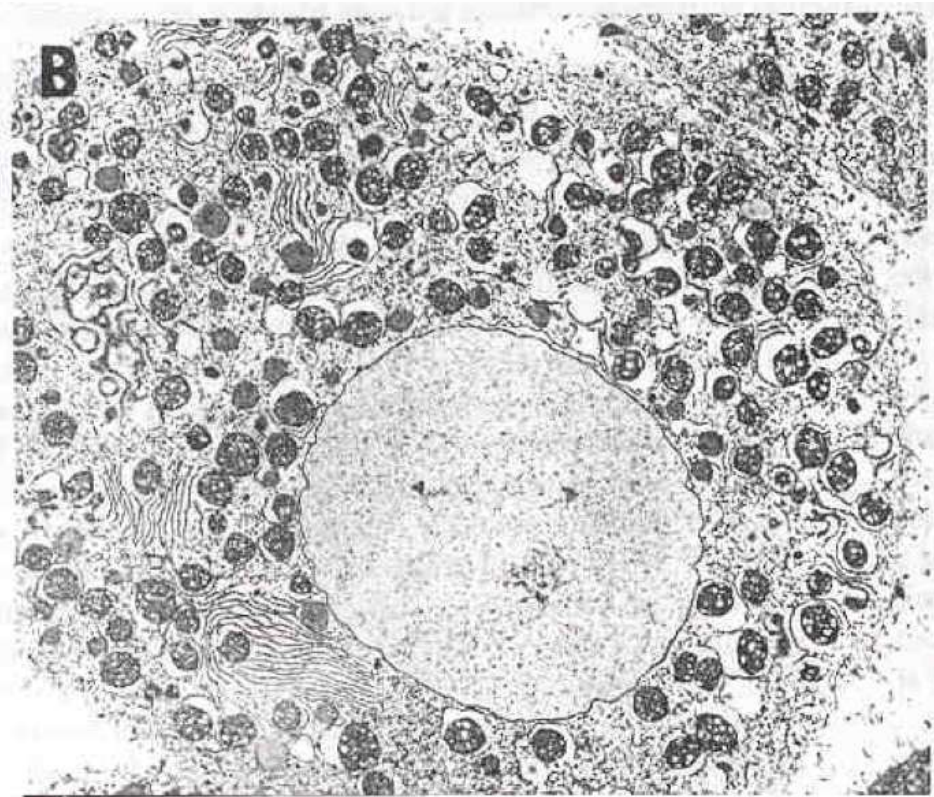


Permeabilized cells provide control over mitochondrial substrate provision

Intact cell



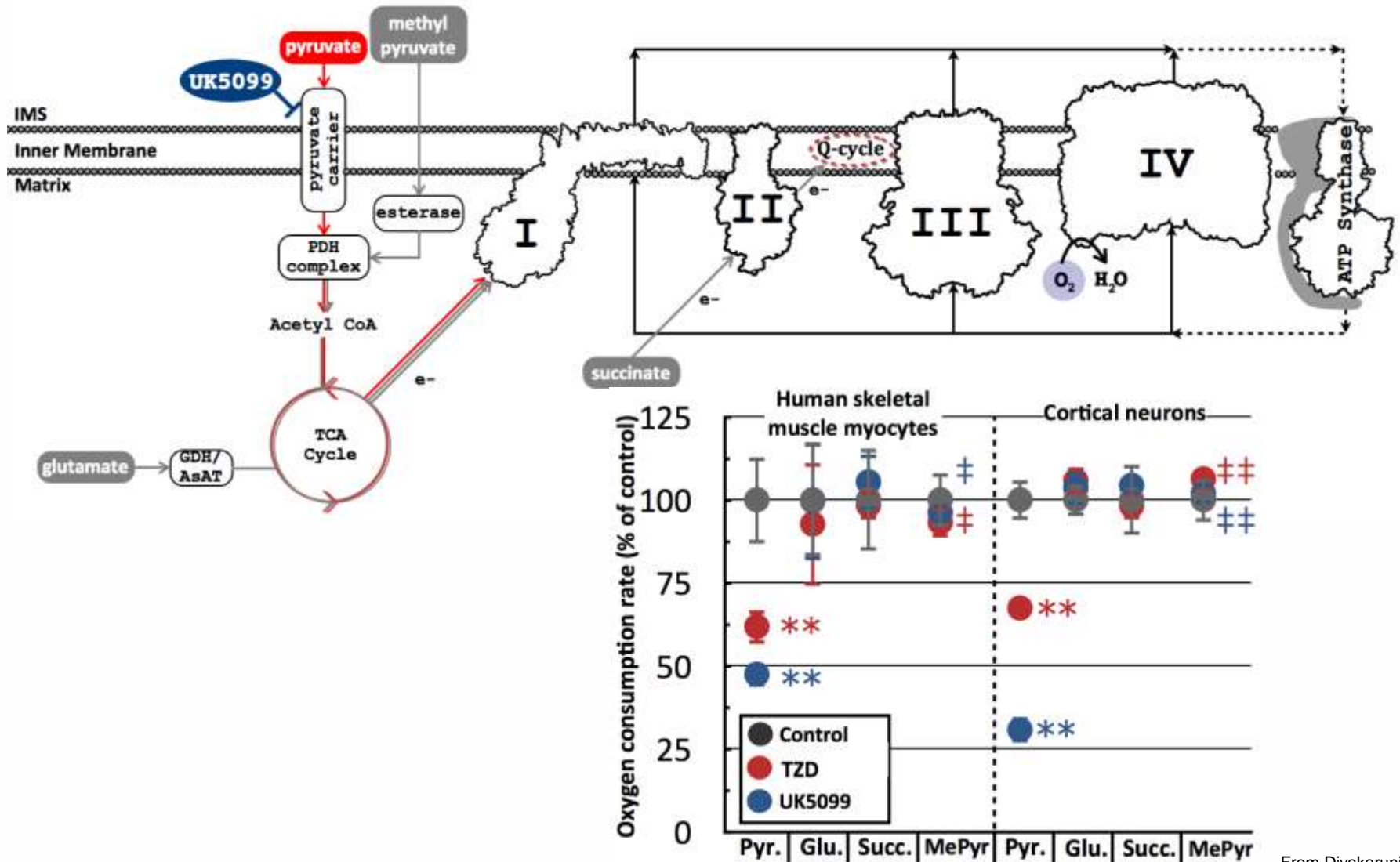
Permeabilized cell



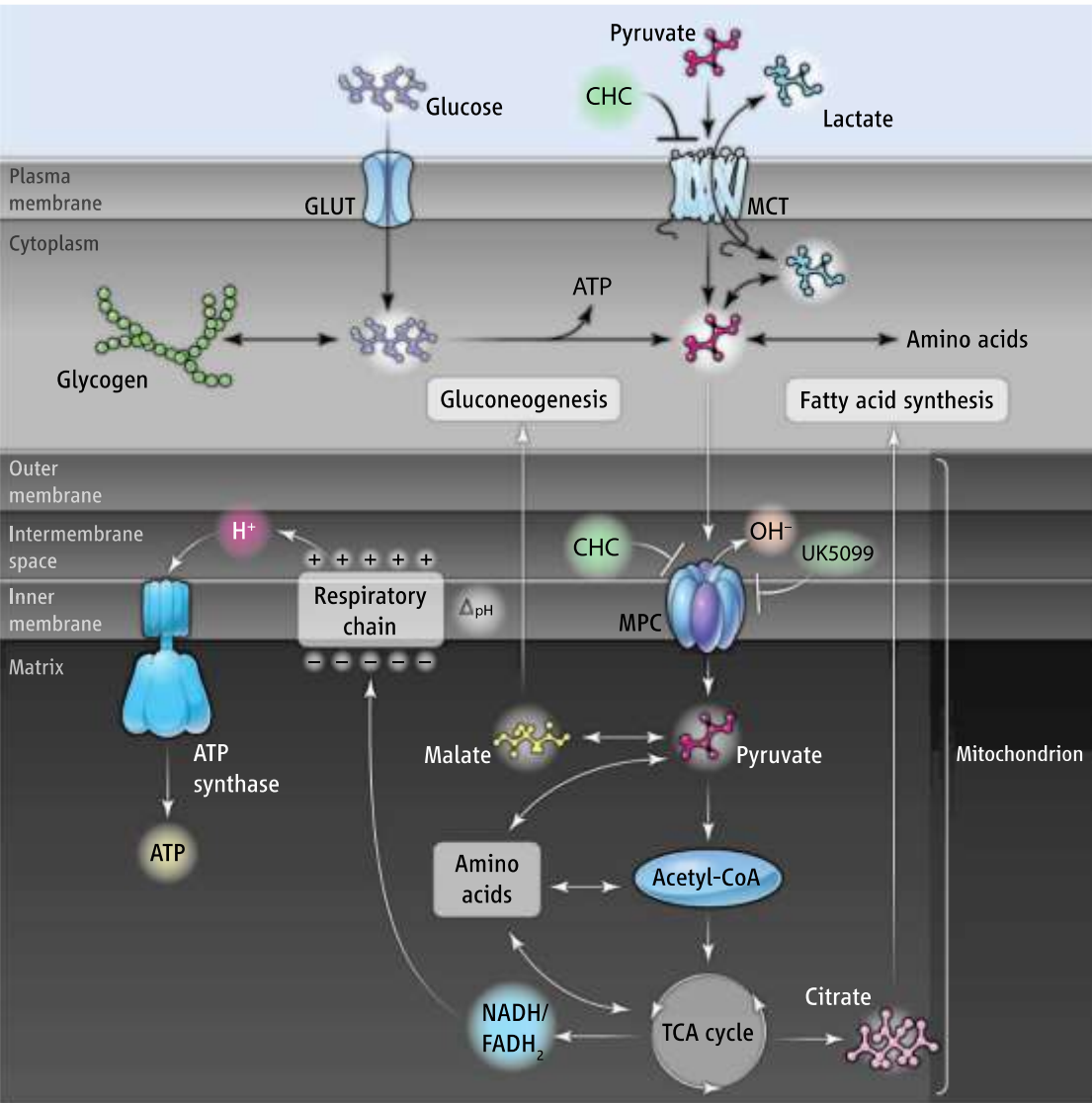
Validation in Divakaruni...and Murphy (2013) *PNAS*;
Protocols in Divakaruni...and Murphy (2014) *Curr. Prot. Tox.*

Image from Fiskum...and Lehninger (1980) *PNAS*

TZDs acutely inhibit mitochondrial pyruvate uptake

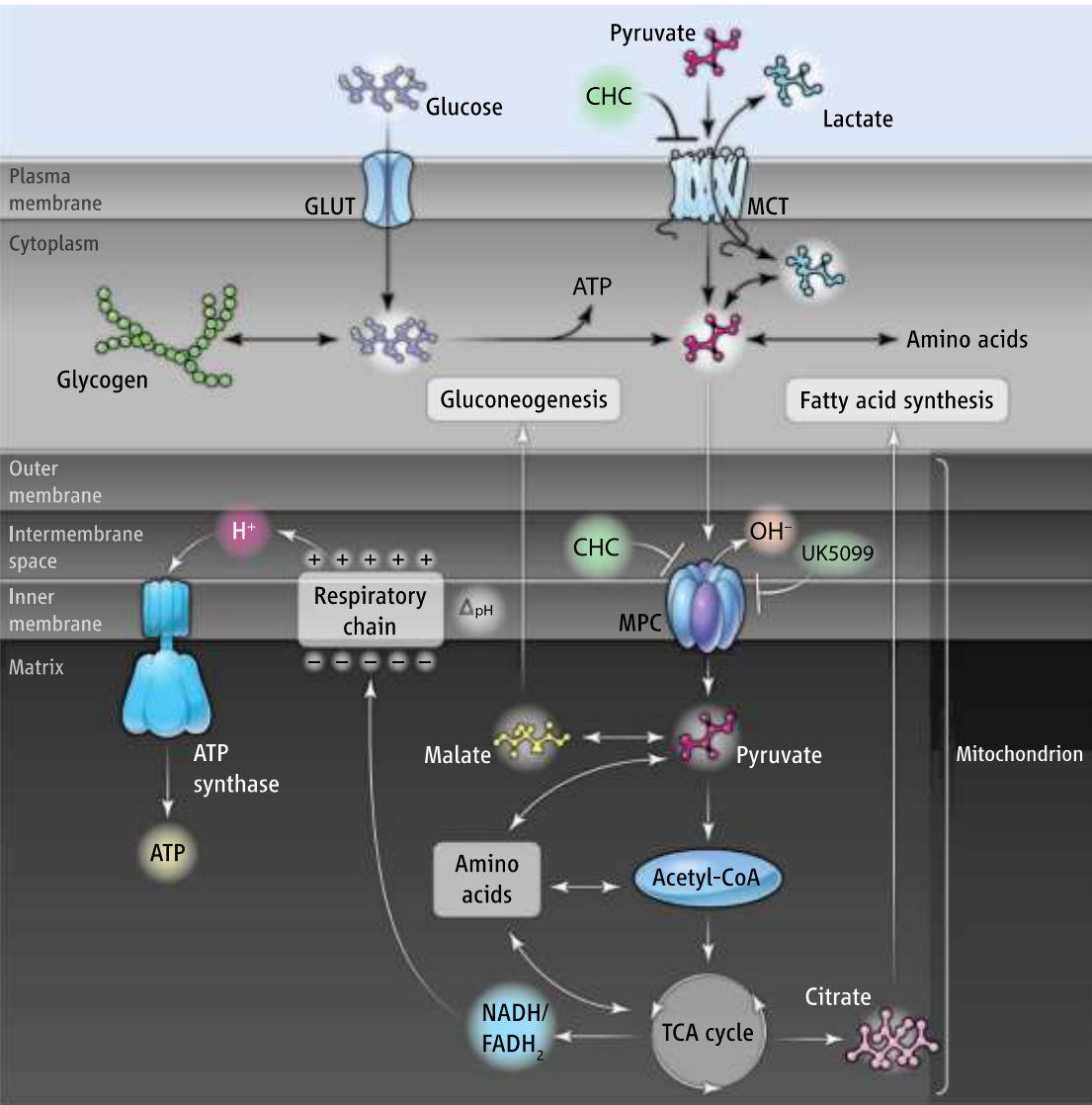


The mitochondrial pyruvate carrier (MPC)



- MPC Knockdown lowers TZD EC_{50}
- UK5099 reproduces plasma membrane glucose uptake and AMPK phosphorylation
Divakaruni and Murphy (2013) *PNAS*
- ¹²⁵I-pioglitazone binds to MPC
Colca and Kletzein (2013) *PLoS ONE*

The mitochondrial pyruvate carrier (MPC)



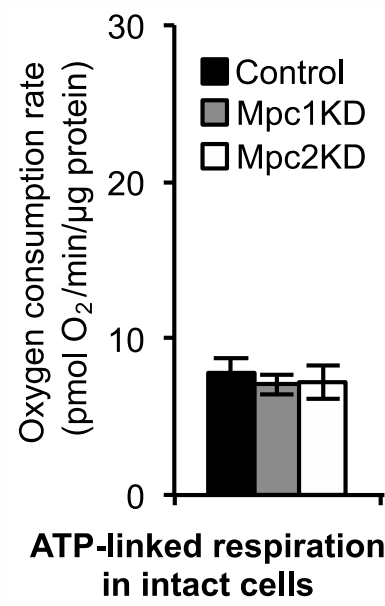
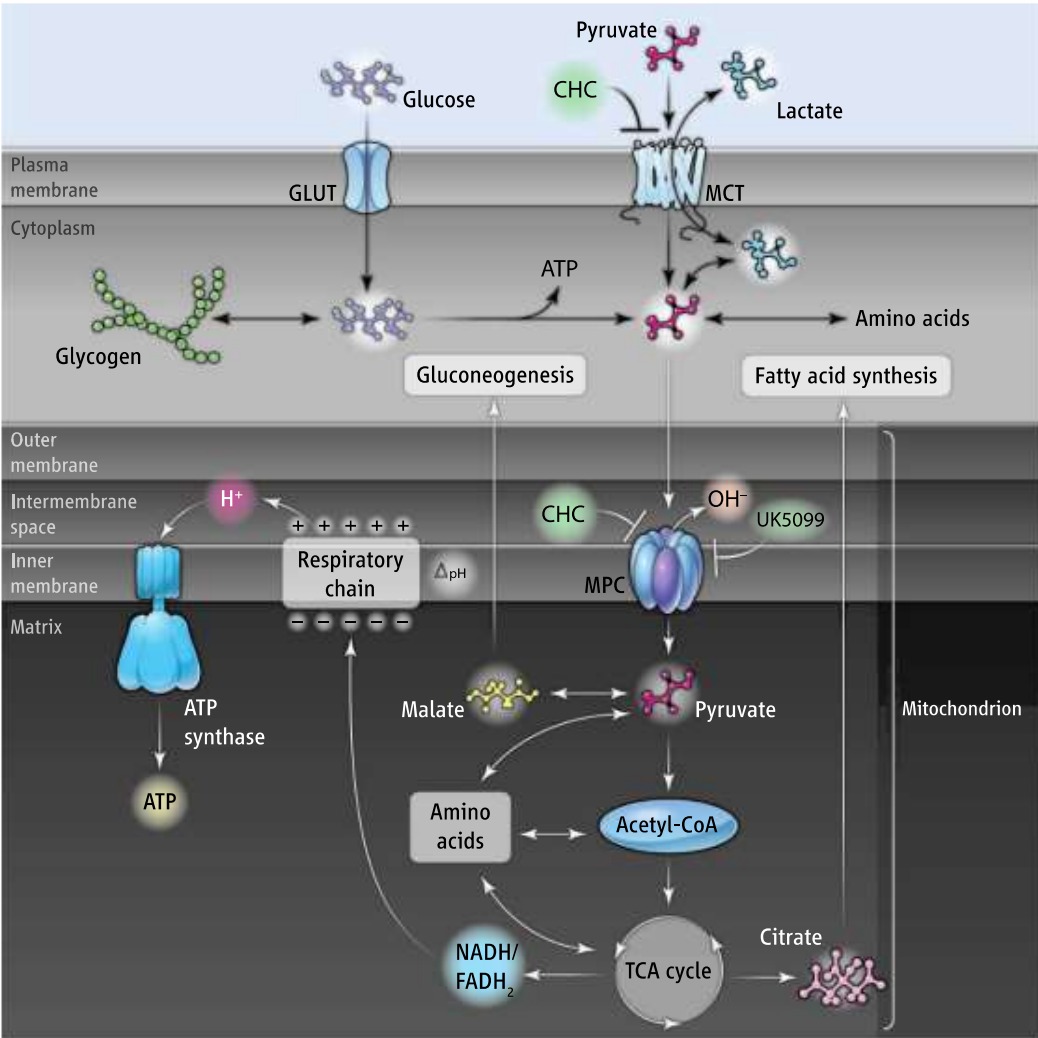
- MPC Knockdown lowers TZD EC₅₀
- UK5099 reproduces plasma membrane glucose uptake and AMPK phosphorylation
Divakaruni and Murphy (2013) *PNAS*
- ¹²⁵I-pioglitazone binds to MPC
Colca and Kletzein (2013) *PLoS ONE*

A METABOLOMICS APPROACH?

- (1) Metabolite changes are not always direct indications of pathways
- (2) Attributing effects to transcriptional agonism vs. MPC inhibition

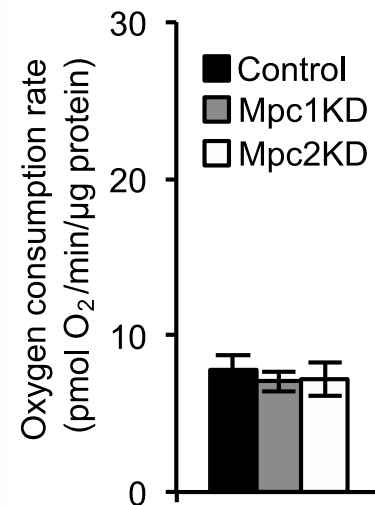
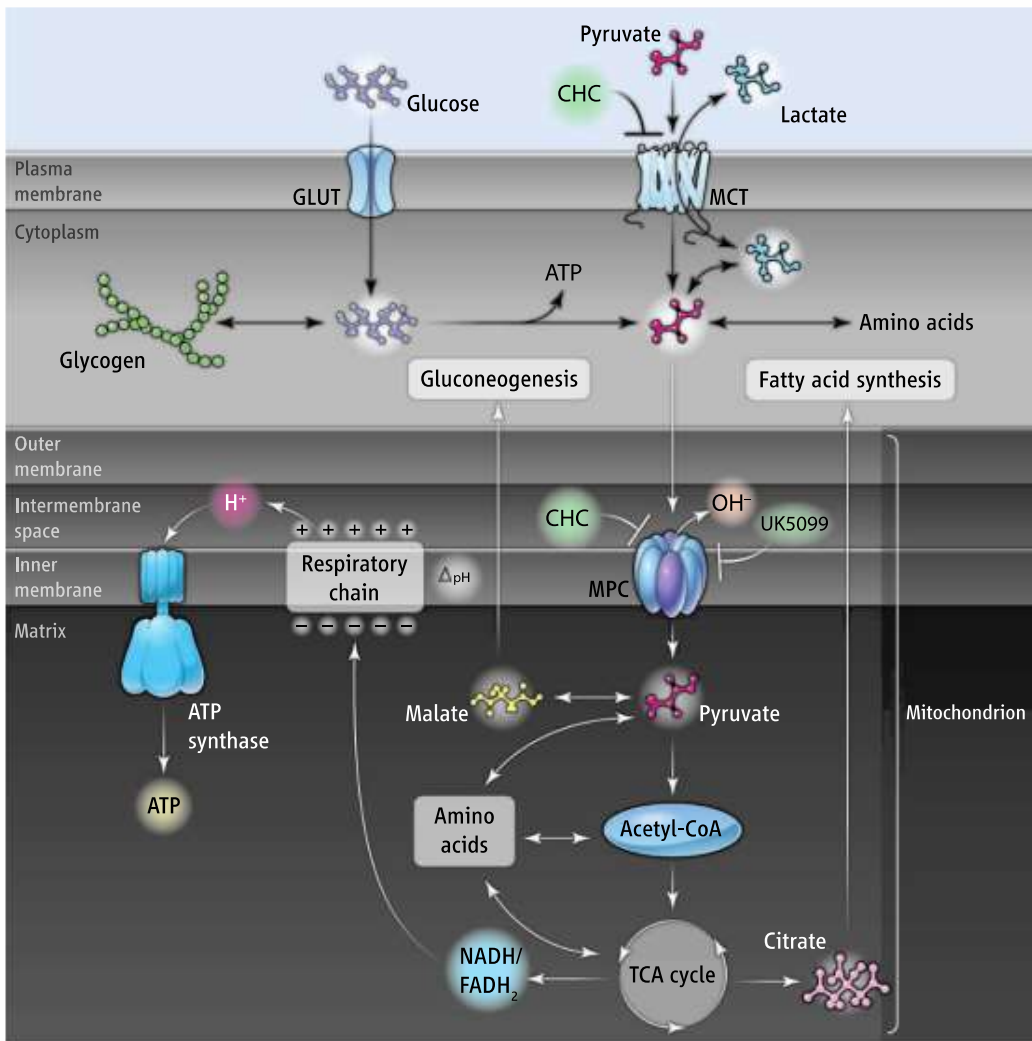
Using stable knockdown to understand the cellular response to reduced MPC activity

**KNOCKING DOWN MPC COMPONENTS
IN MYOCYTES DOESN'T ALTER
MITOCHONDRIAL ATP PRODUCTION (!)**

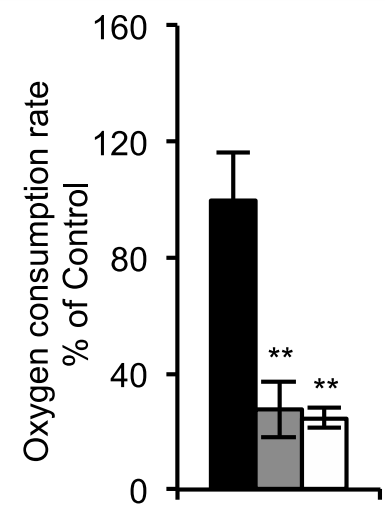


Using stable knockdown to understand the cellular response to reduced MPC activity

**KNOCKING DOWN MPC COMPONENTS
IN MYOCYTES DOESN'T ALTER
MITOCHONDRIAL ATP PRODUCTION (!)**



ATP-linked respiration in intact cells



Pyruvate-driven respiration in permeabilized cells

Metabolism, two ways

**(1) MEASURE RATES OF
INPUTS AND OUTPUTS**

**(2) TAKE A DETAILED
SNAPSHOT IN TIME**

*(Metabolomics, stable isotope
tracing)*

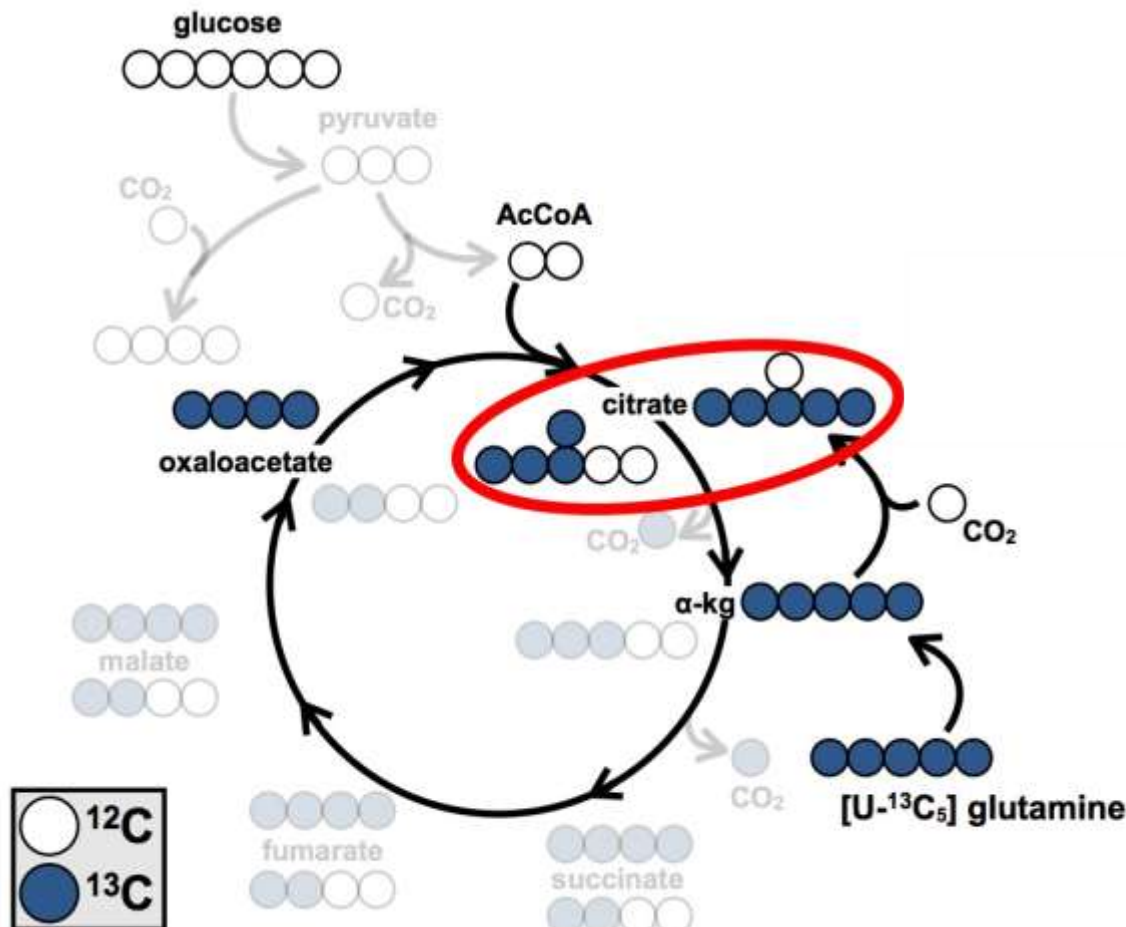
Granular detail from
metabolite abundances

Relative enrichment of
intermediates from labeled
substrates

Combining respiration and stable isotope tracing to fully reveal cell metabolism

Isotopically labeled substrates (i.e. glucose, glutamine, etc.) are offered as tracers

^{13}C labeling of metabolite pools can map the pattern of carbon flux

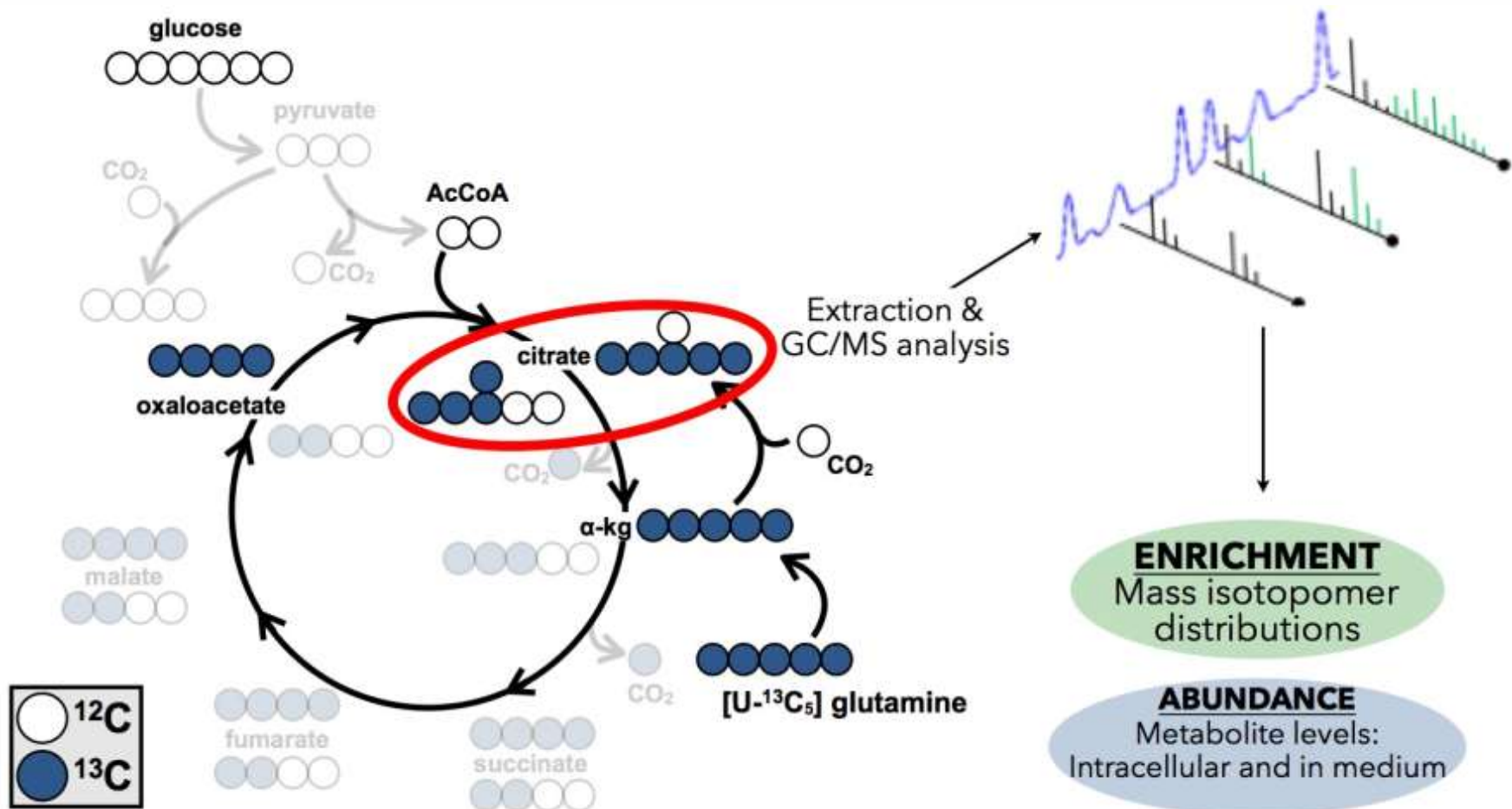


For Research Use Only. Not for use in diagnostic procedures.

Combining respiration and stable isotope tracing to fully reveal cell metabolism

Isotopically labeled substrates (i.e. glucose, glutamine, etc.) are offered as tracers

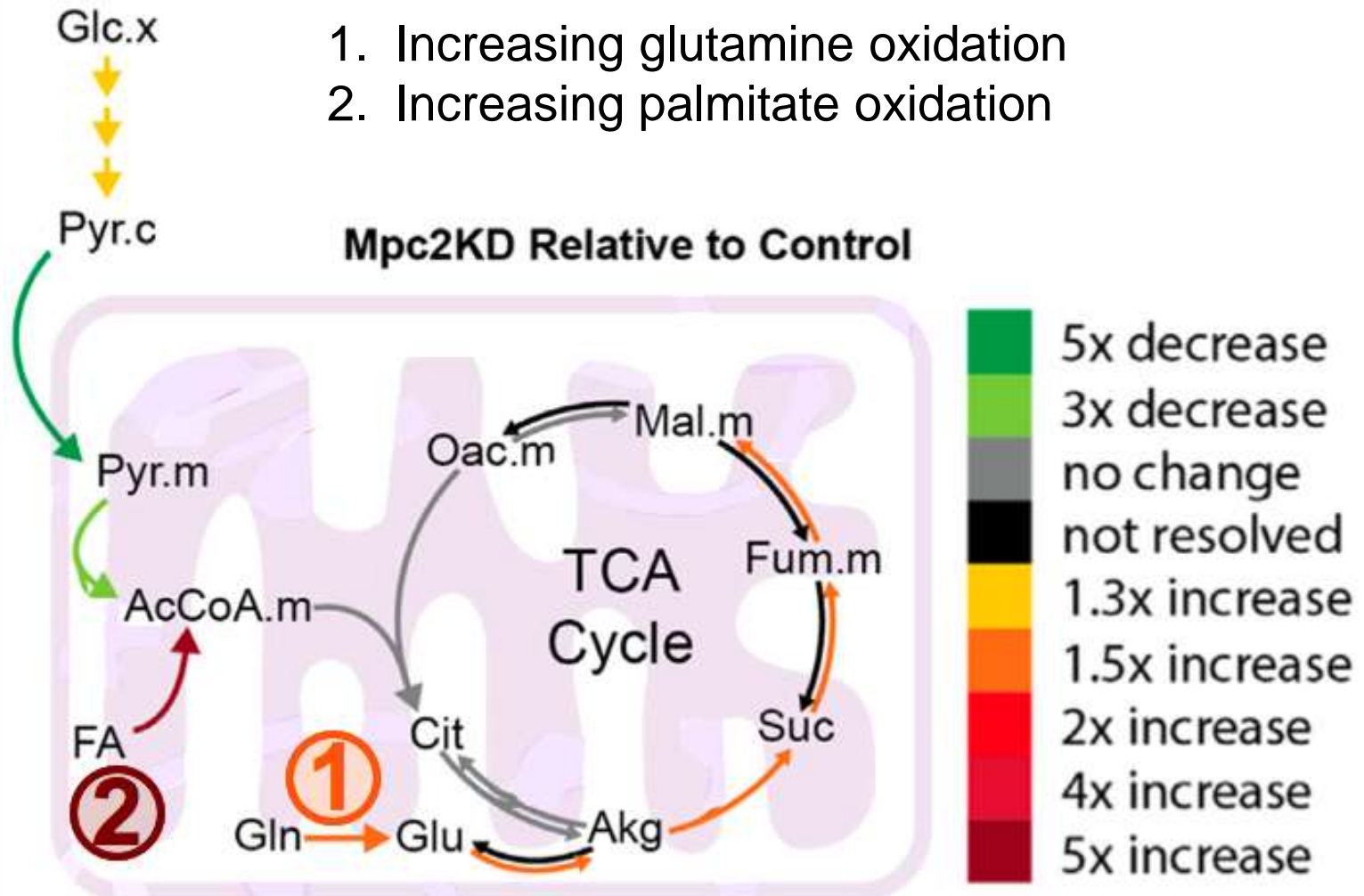
^{13}C labeling of metabolite pools can map the pattern of carbon flux



The MPC is a switch that can potentiate the oxidation of amino acids and fatty acids

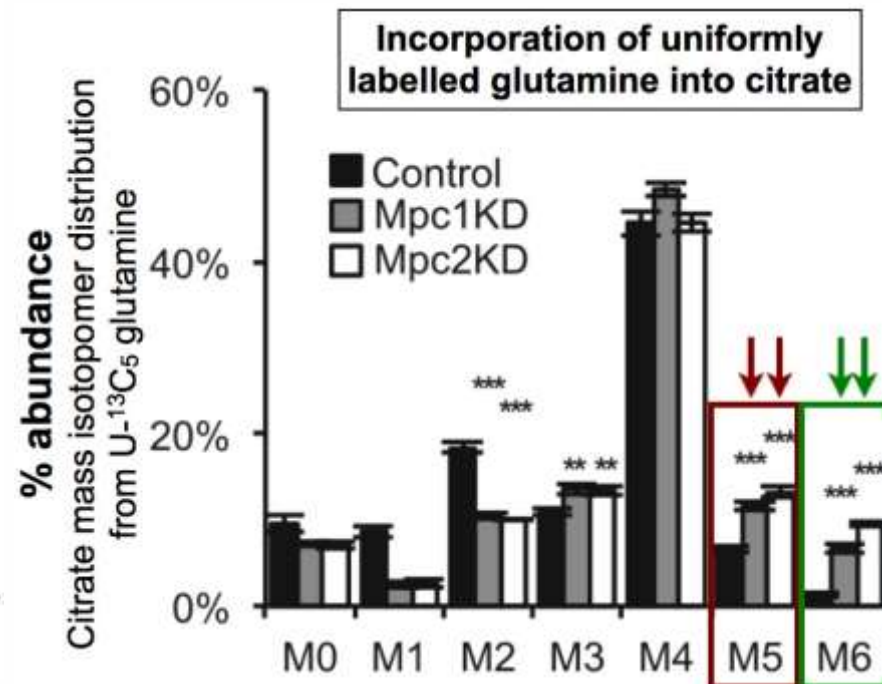
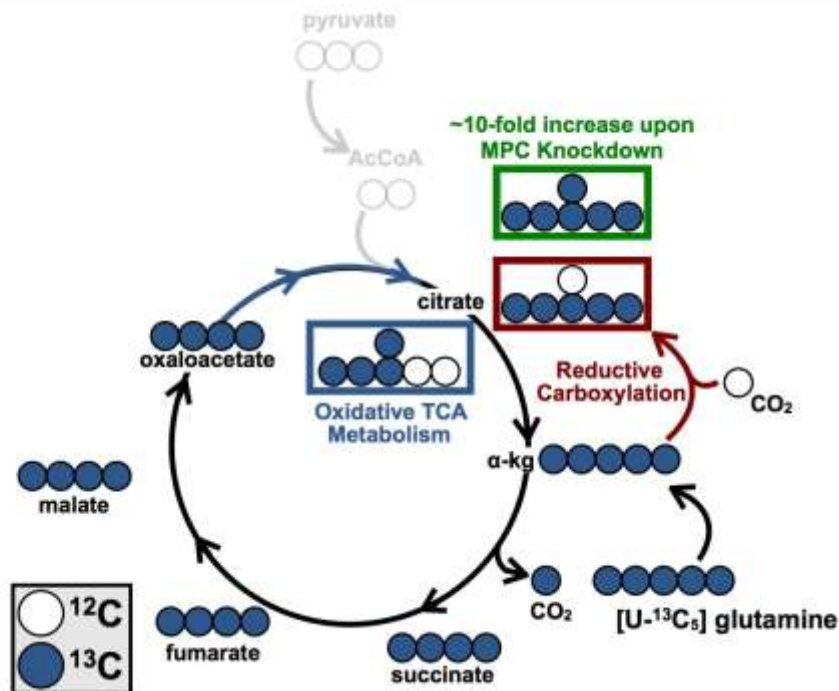
MYOCYTES ADAPT TO DECREASED MITOCHONDRIAL PYRUVATE UPTAKE BY:

1. Increasing glutamine oxidation
2. Increasing palmitate oxidation



Stable isotope tracing can reveal information that respirometry cannot

For research use only. Not for use in diagnostic procedures.



For Research Use Only. Not for use in diagnostic procedures.

Malic enzyme activity increases upon MPC knockdown

Glutamine Oxidation Maintains the TCA Cycle and Cell Survival during Impaired Mitochondrial Pyruvate Transport

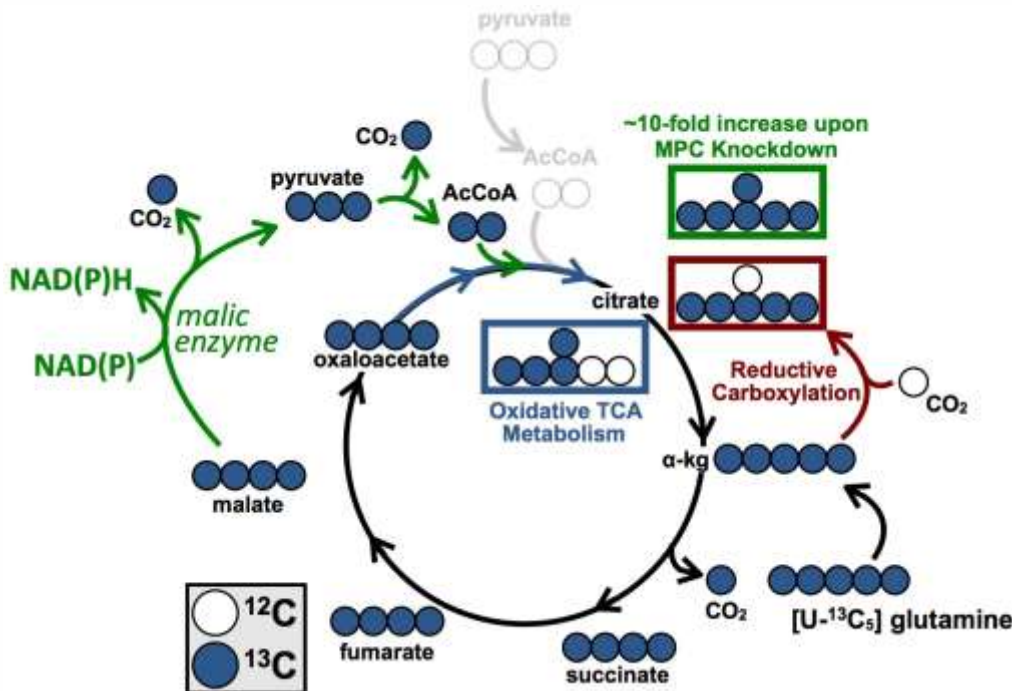
Chendong Yang,¹ Bookyung Ko,¹ Christopher T. Hensley,¹ Lei Jiang,¹ Aja T. Wasti,^{1,2} Jiyeon Kim,¹ Jessica Sudderth,¹ Maria Antonietta Calvaruso,¹ Lloyd Lumata,² Matthew Mitsche,² Jared Rutter,² Matthew E. Merritt,³ and Ralph J. DeBerardinis^{1,2,4,*}

Molecular Cell 56 November 6, 2014

Regulation of Substrate Utilization by the Mitochondrial Pyruvate Carrier

Nathaniel M. Vacanti,¹ Ajit S. Divakaran,² Courtney R. Green,¹ Seth J. Parker,¹ Robert R. Henry,^{3,4} Theodore P. Claraldi,^{3,4} Anne N. Murphy,² and Christian M. Metallo^{1,2}

Molecular Cell 56 November 6, 2014



Citrate mass isotopomer distribution from U-¹³C₅ glutamine

Incorporation of uniformly labelled glutamine into citrate

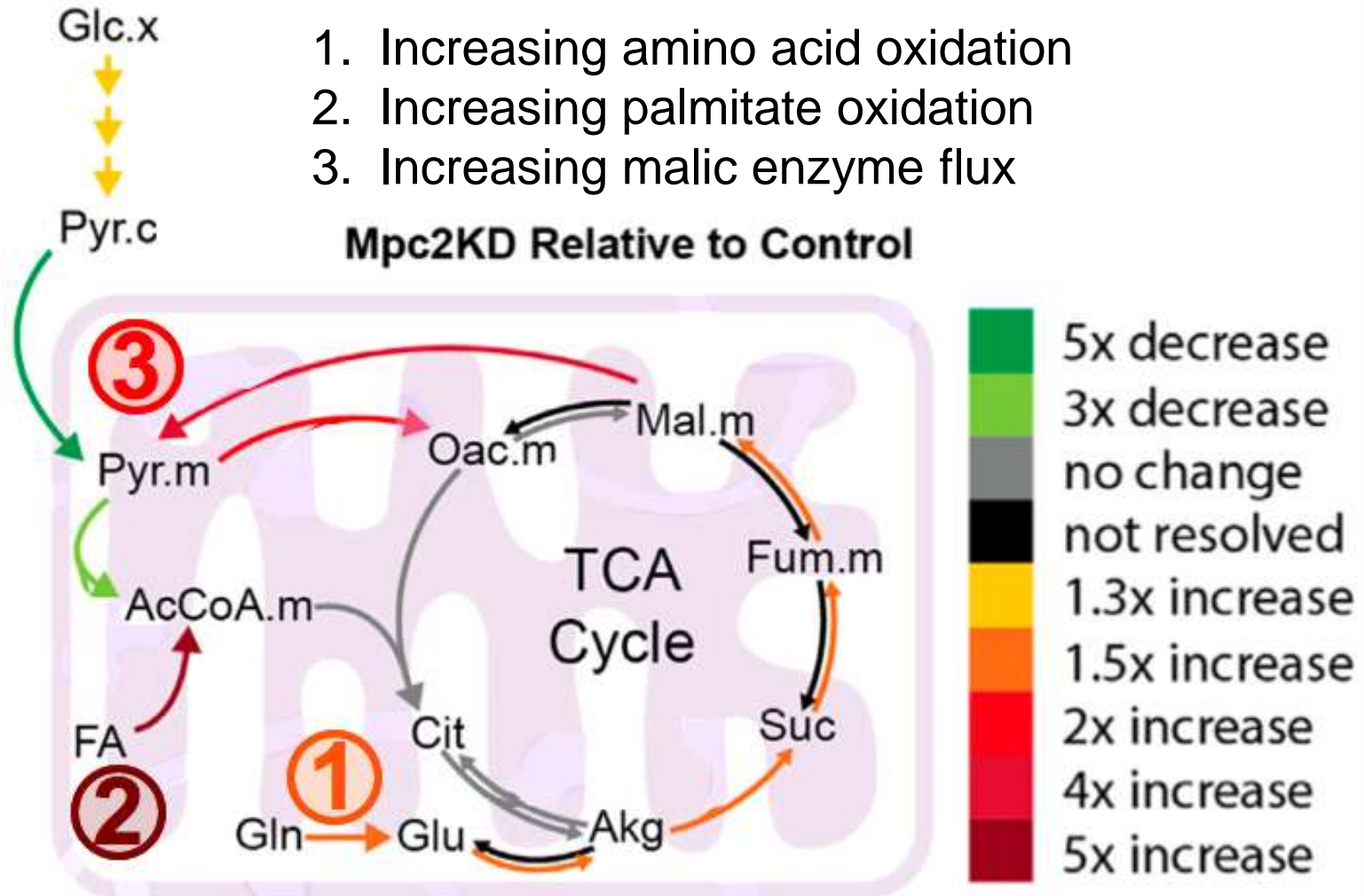
Control
Mpc1KD
Mpc2KD

M0 M1 M2 M3 M4 M5 M6

The MPC is a “switch” that can potentiate metabolic flexibility

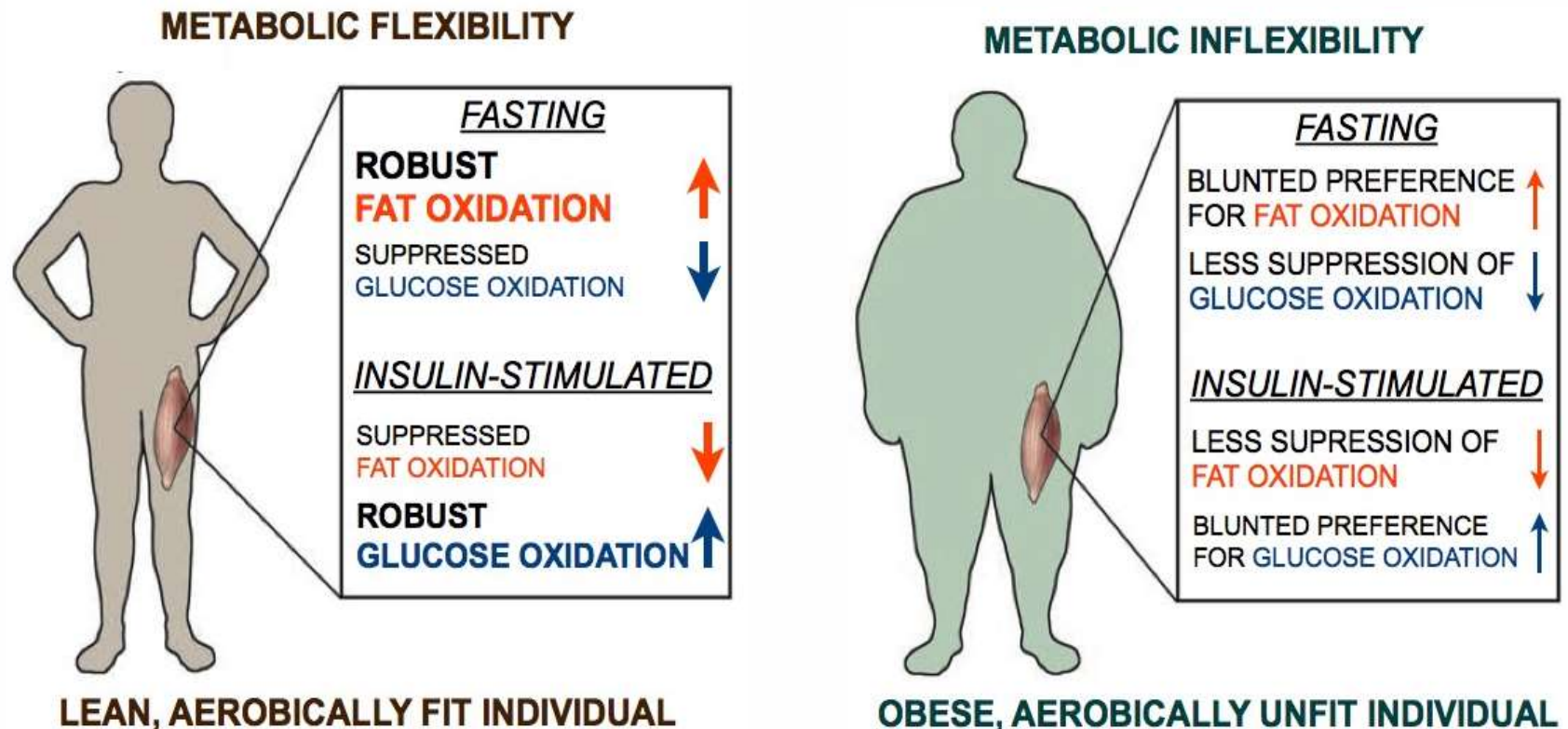
MYOCYTES ADAPT TO DECREASED MITOCHONDRIAL PYRUVATE UPTAKE BY:

1. Increasing amino acid oxidation
2. Increasing palmitate oxidation
3. Increasing malic enzyme flux



The MPC is a “switch” that can potentiate metabolic flexibility

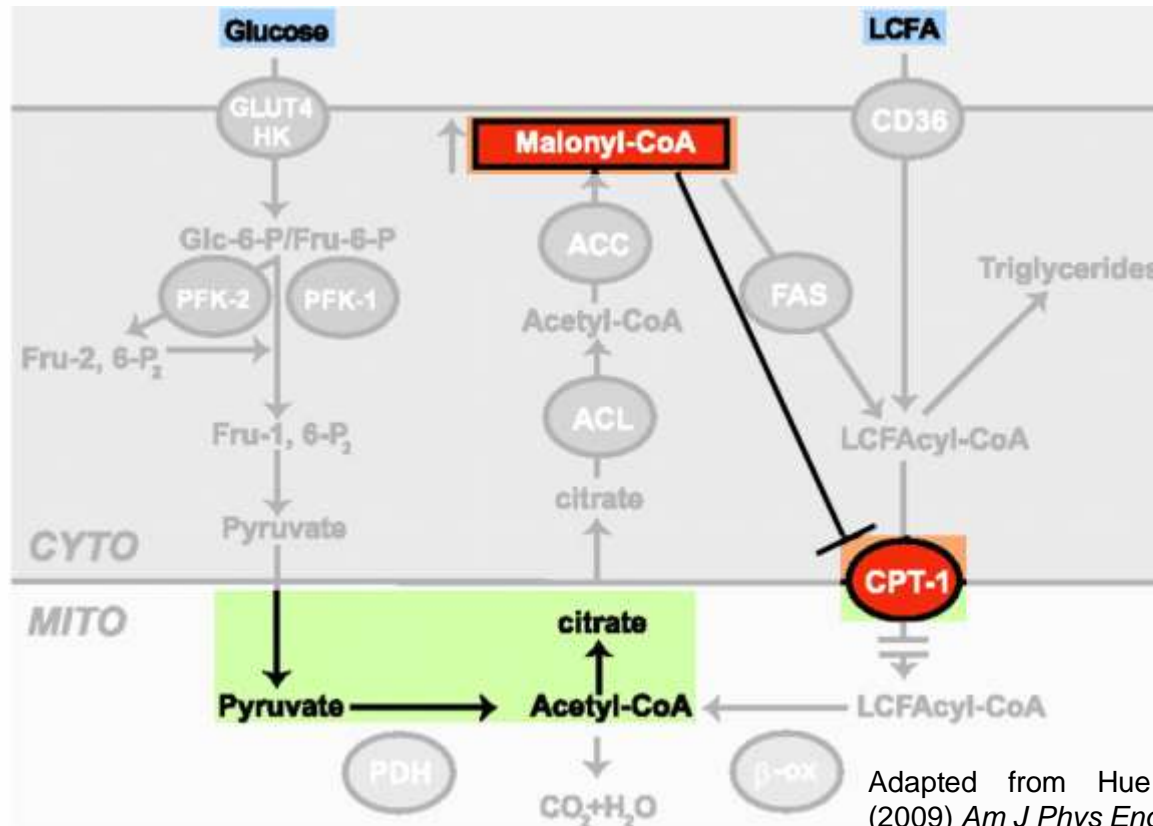
Metabolic flexibility is a characteristic of healthy skeletal muscle



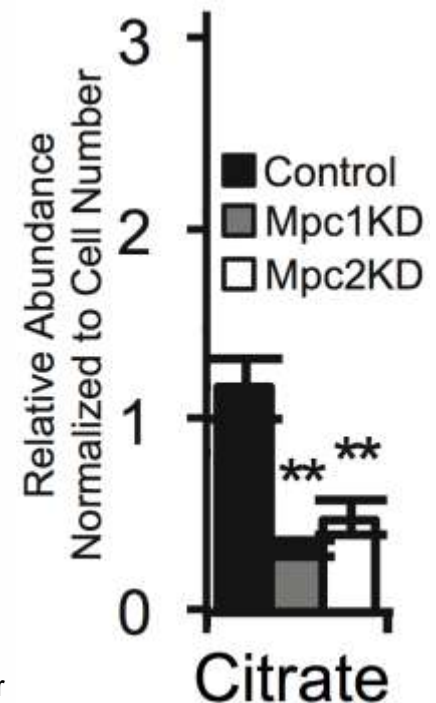
The MPC is a “switch” that can potentiate metabolic flexibility

MPC inhibition potentiates fatty acid oxidation (by reducing malonyl CoA?)

An inability to increase lipid oxidation when faced with high supply can cause increased muscle fat accumulation and insulin resistance



Adapted from Hue and Taegtmeyer (2009) *Am J Phys End Metab*



From Vacanti,...and Metallo (2014) *Mol. Cell*

Is there a metabolic component to neurodegenerative disease?

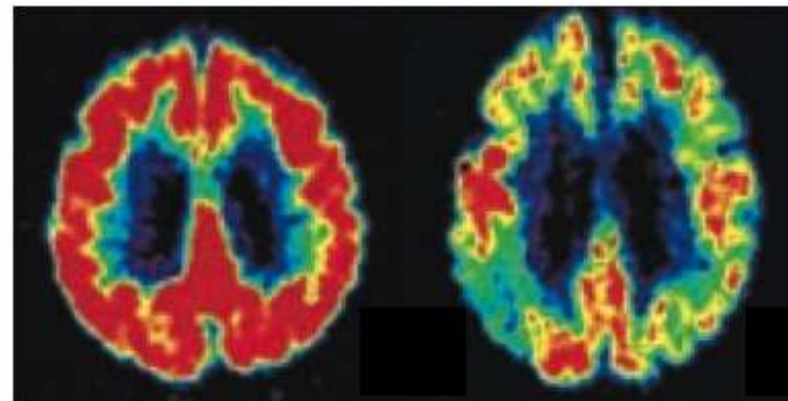
TYPE 2 DIABETES IS A STRONG RISK FACTOR FOR AD

SHARED PATHOLOGY BETWEEN T2D AND AD

- Insulin resistance
- Amyloidosis
- Oxidative stress
- Cognitive impairment
- Neural tissue atrophy

Healthy brain

Mild-to-moderate AD



Is there a metabolic component to neurodegenerative disease?

TYPE 2 DIABETES IS A STRONG RISK FACTOR FOR AD

SHARED PATHOLOGY BETWEEN T2D AND AD

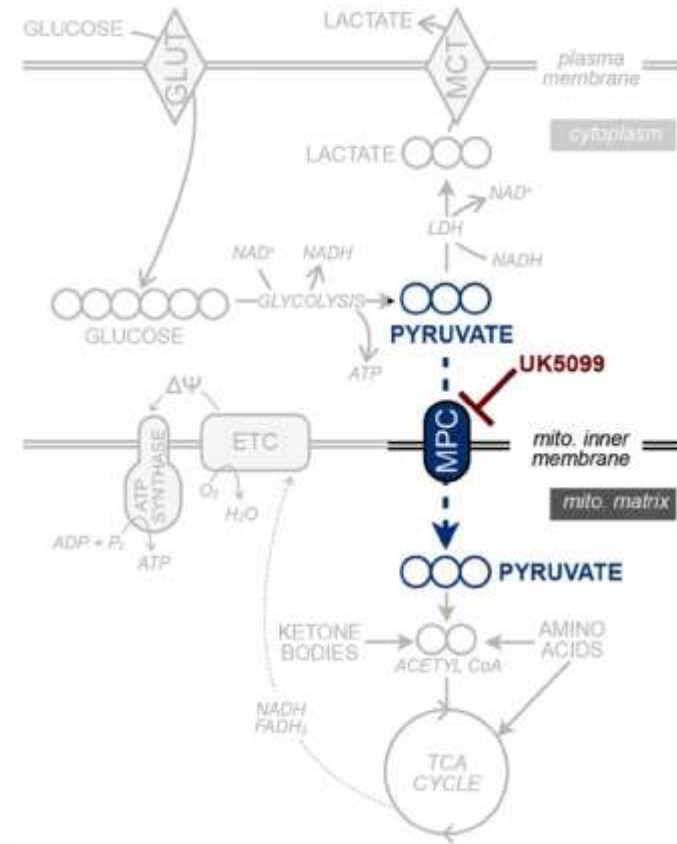
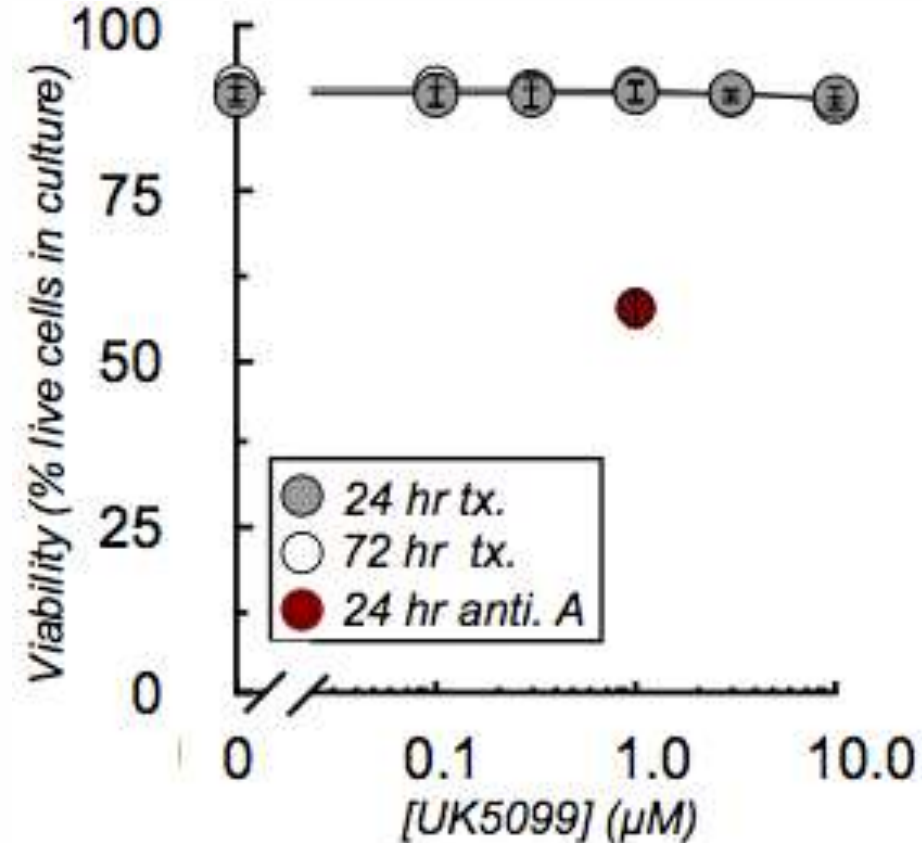
- Insulin resistance
- Amyloidosis
- Oxidative stress
- Cognitive impairment
- Neural tissue atrophy

REPURPOSING OF ANTI-DIABETIC THERAPIES FOR AD

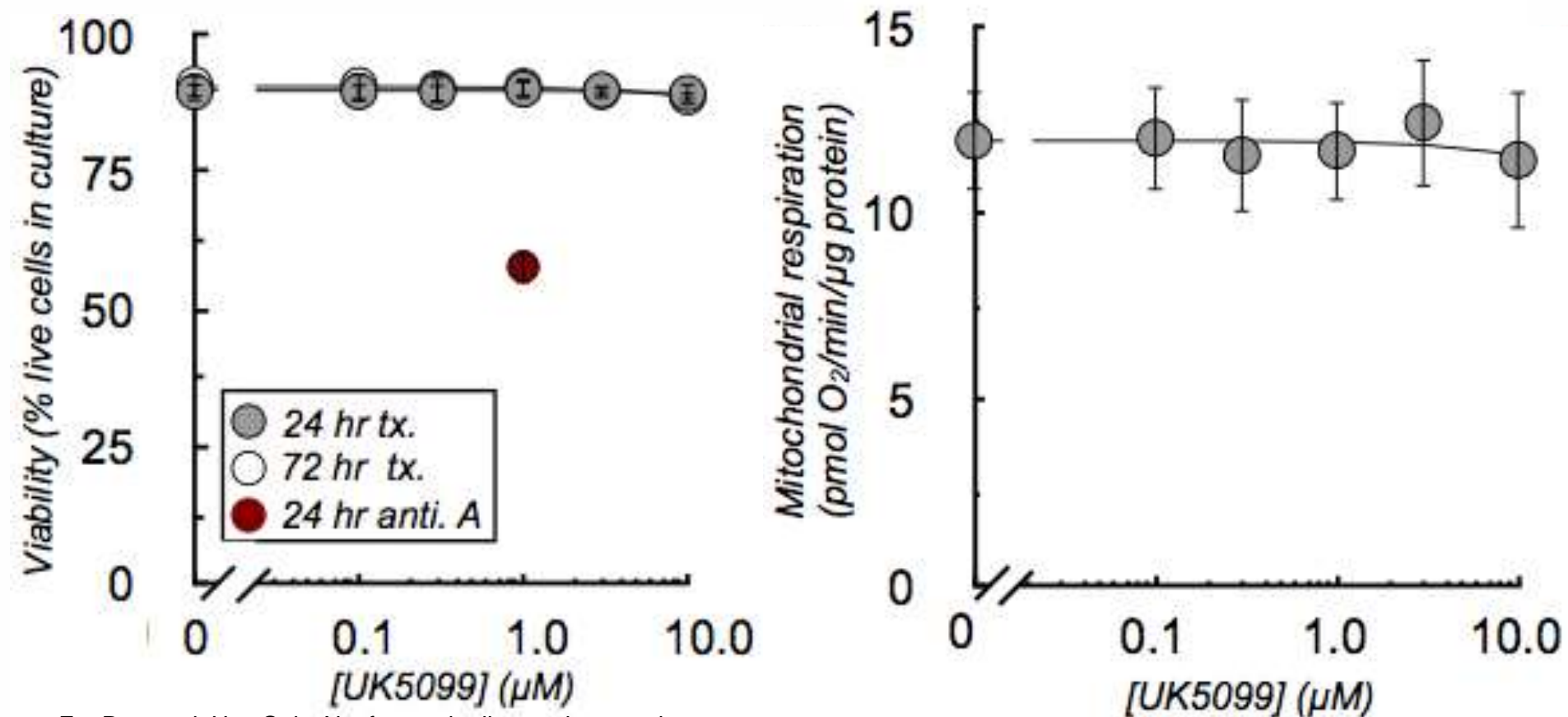
- Intranasal insulin delivery
- Incretin analogs
- Thiazolidinediones (TZDs)



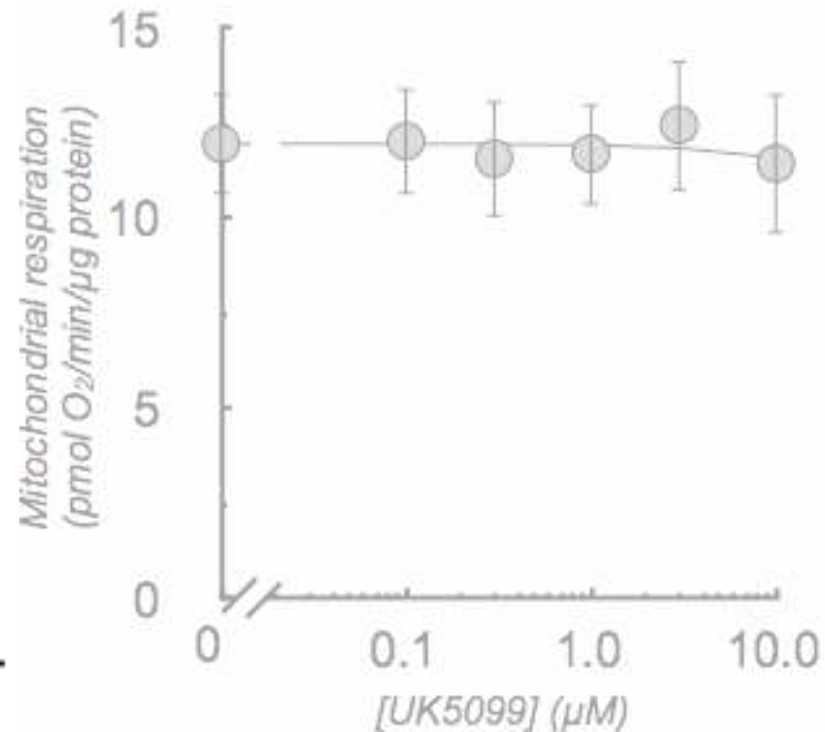
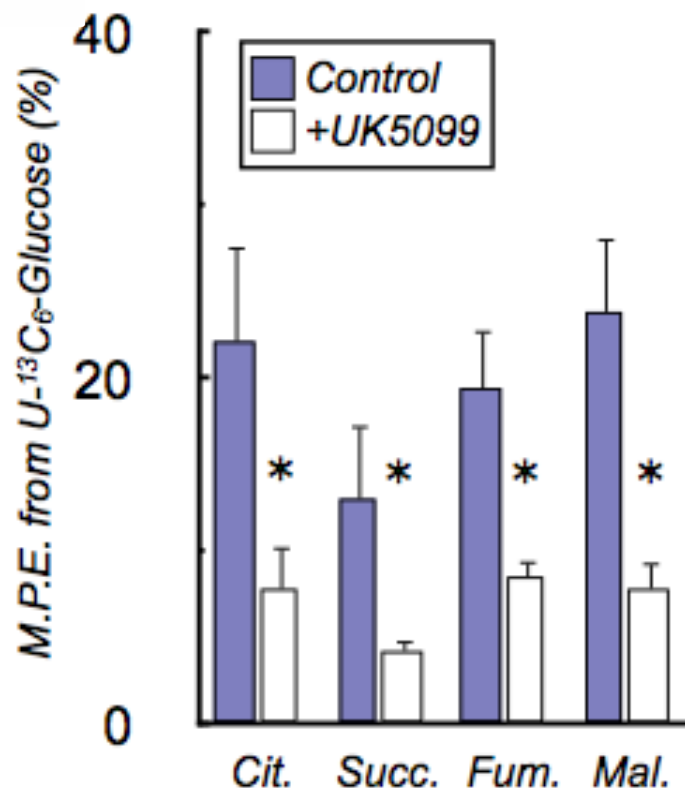
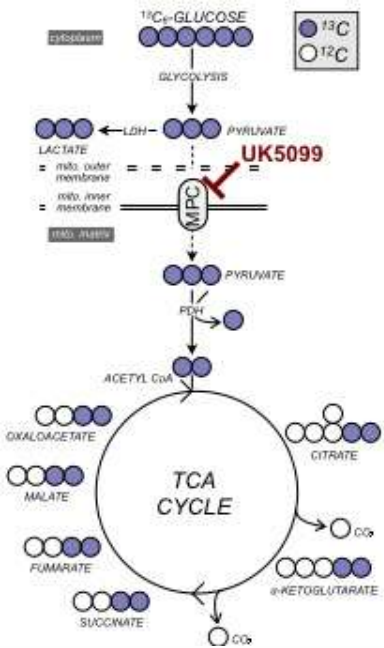
Viability is unchanged despite reduced MPC activity



Global bioenergetics unchanged despite reduced MPC activity

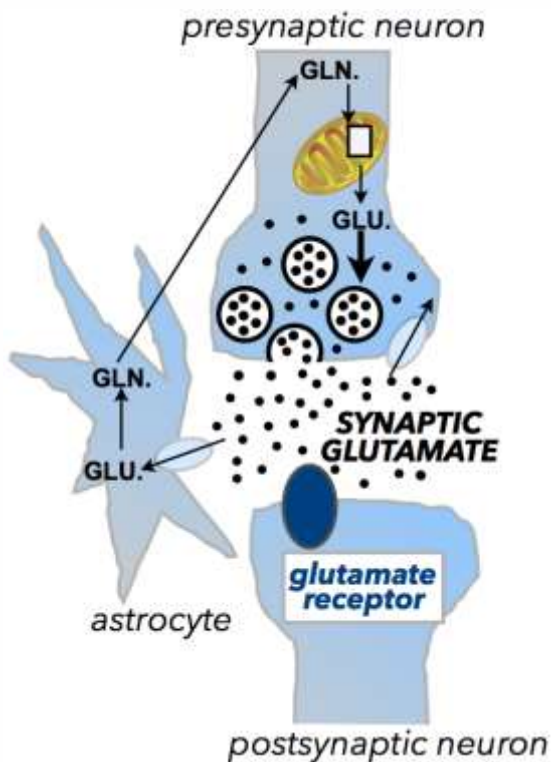


Global bioenergetics unchanged despite reduced MPC activity



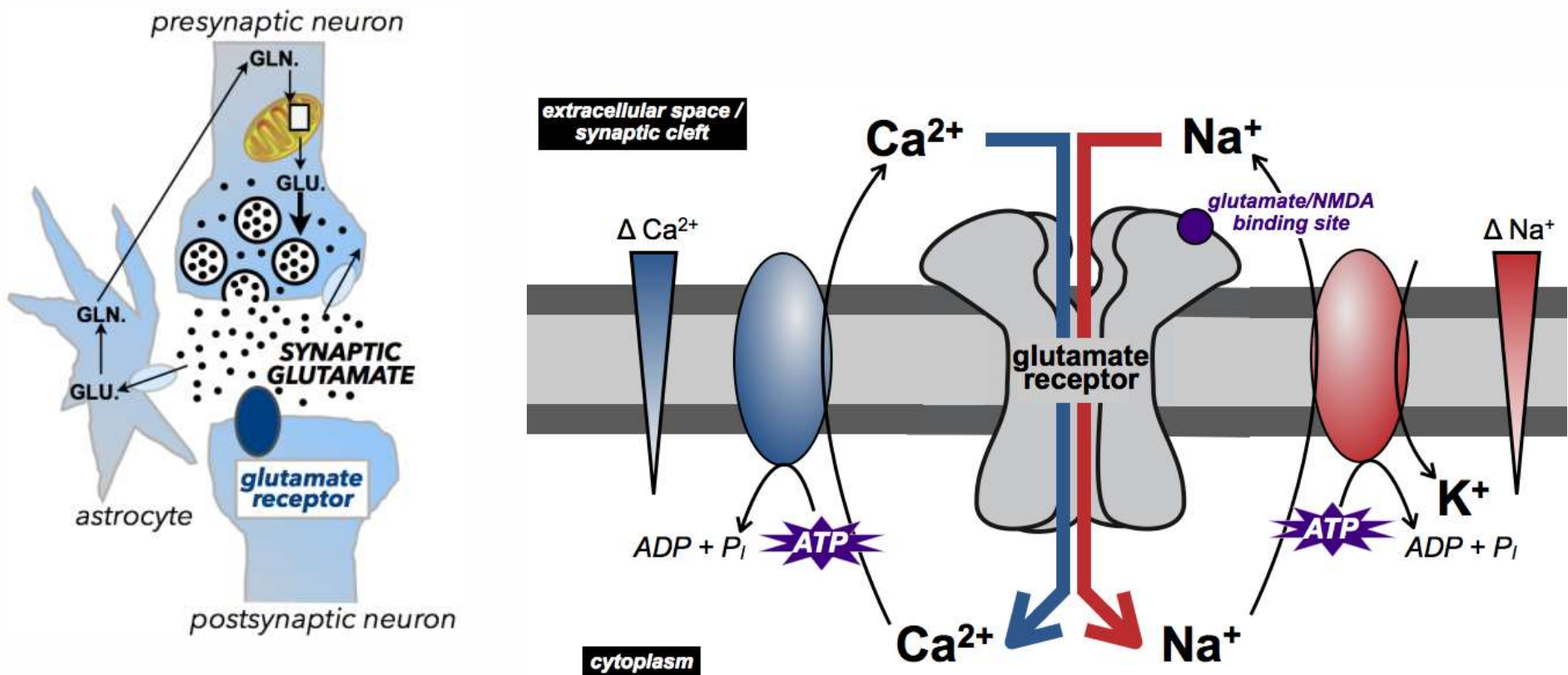
Challenging neurons by stimulating glutamate receptors

Glutamate is the dominant excitatory neurotransmitter in the brain
(balanced by inhibitory GABA)



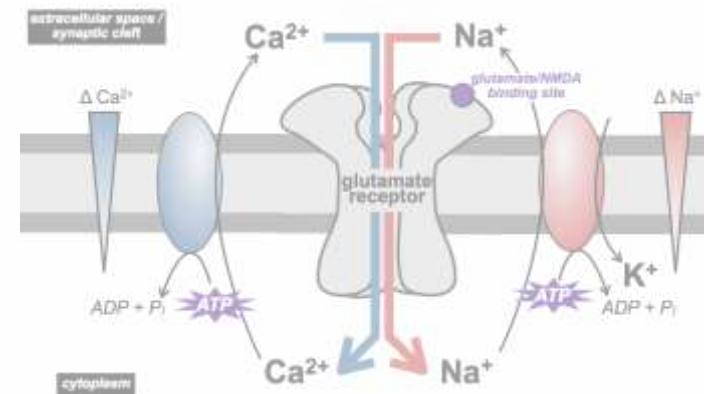
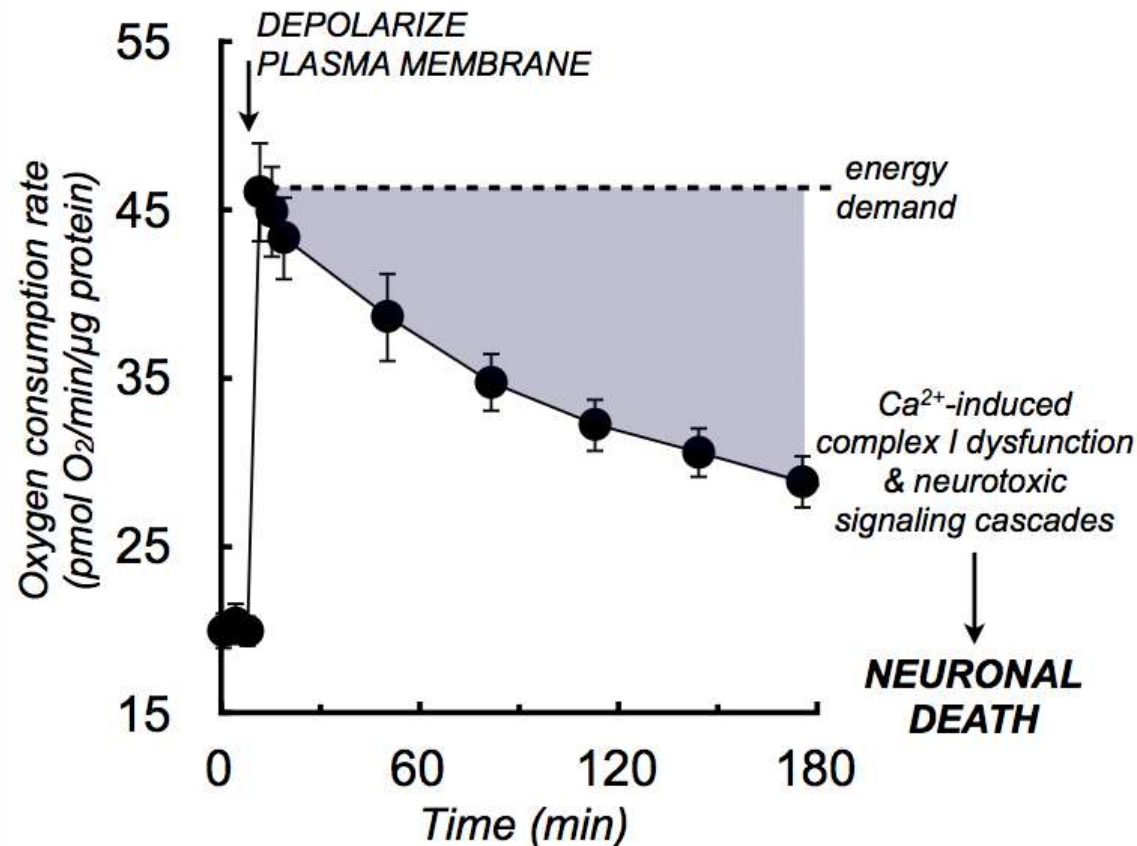
Challenging neurons by stimulating glutamate receptors

Glutamatergic neurotransmission comes at a tremendous energetic cost



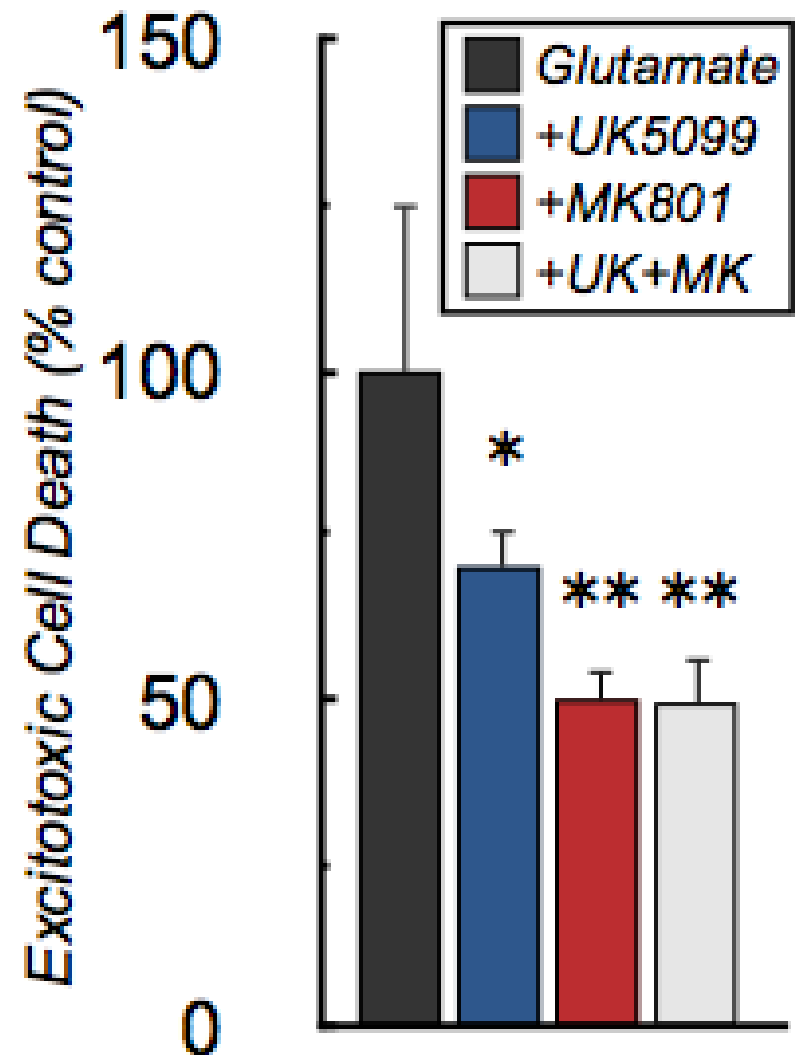
Challenging neurons by stimulating glutamate receptors

Under conditions of energetic stress or de-energization (stroke, epilepsy, some chronic neurodegeneration), extracellular glutamate rises to toxic levels



MPC inhibition protects from excitotoxic cell death (!)

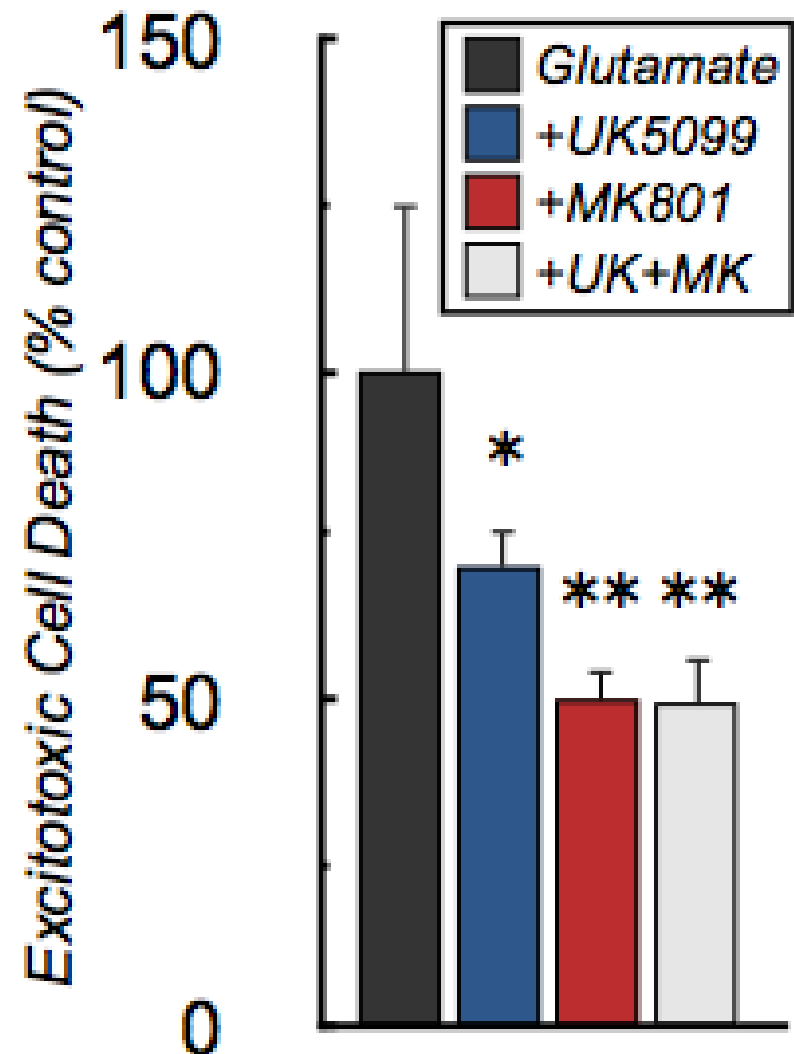
In rich culture medium, inhibiting mitochondrial pyruvate metabolism in neurons protects from excitotoxic neuronal death



MPC inhibition protects from excitotoxic cell death (!)

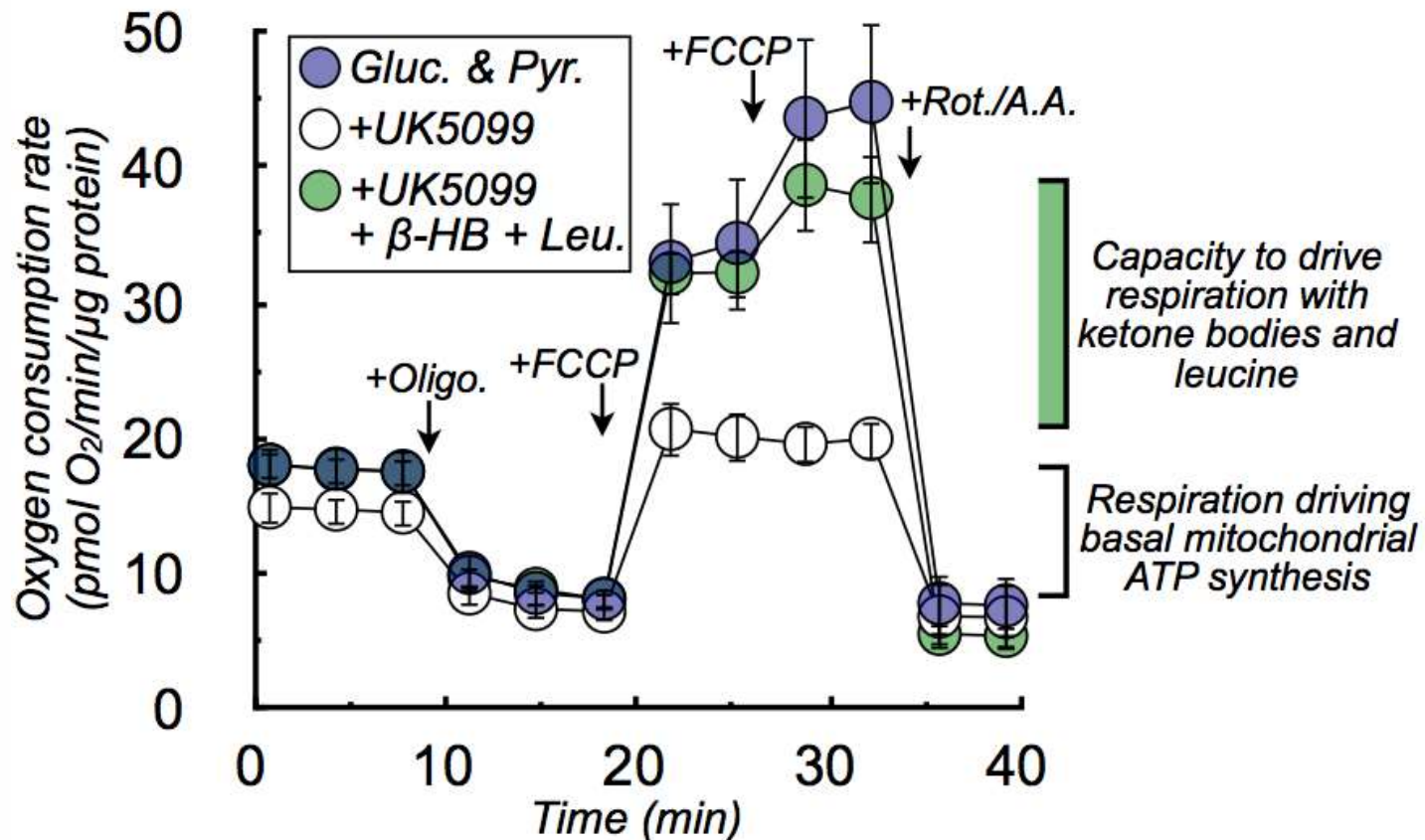
In rich culture medium, inhibiting mitochondrial pyruvate metabolism in neurons protects from excitotoxic neuronal death

How has MPC inhibition rewired neuronal metabolism in a way that is neuroprotective?

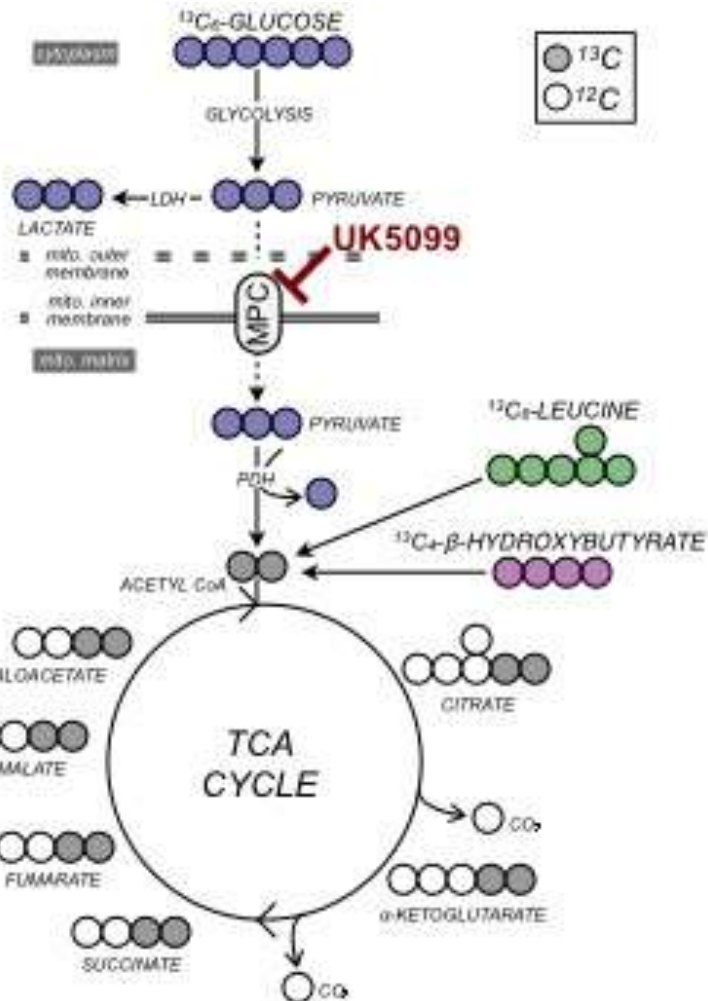


Bioenergetic analysis shows neurons have a huge capacity to oxidize ketone bodies and BCAAs...

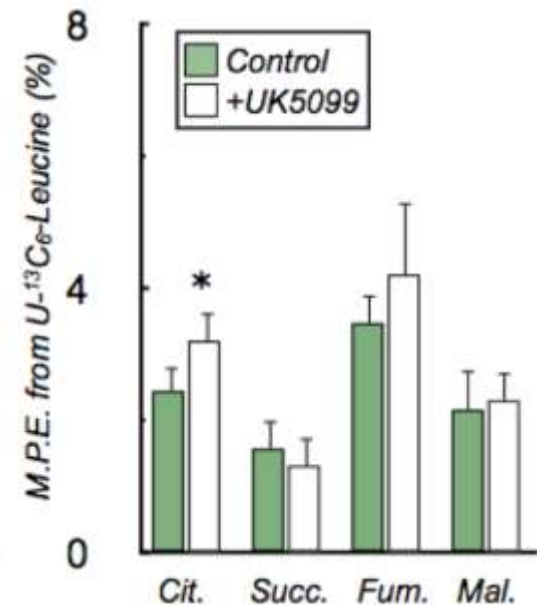
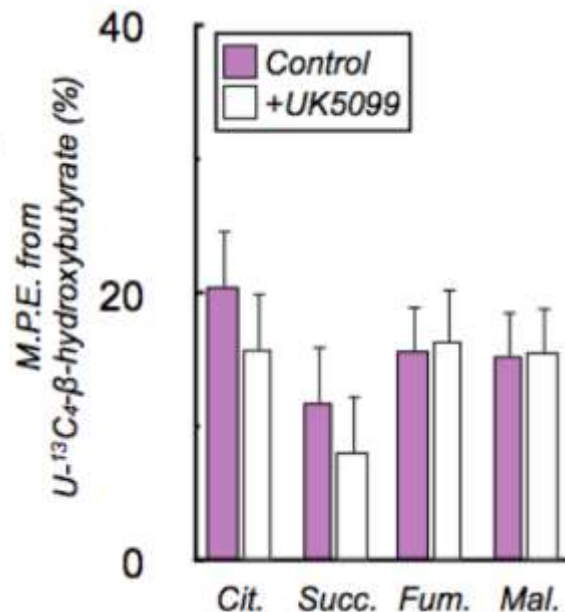
Acute challenge in simple salts medium where substrate preference is assumed *a priori* by the experimenter



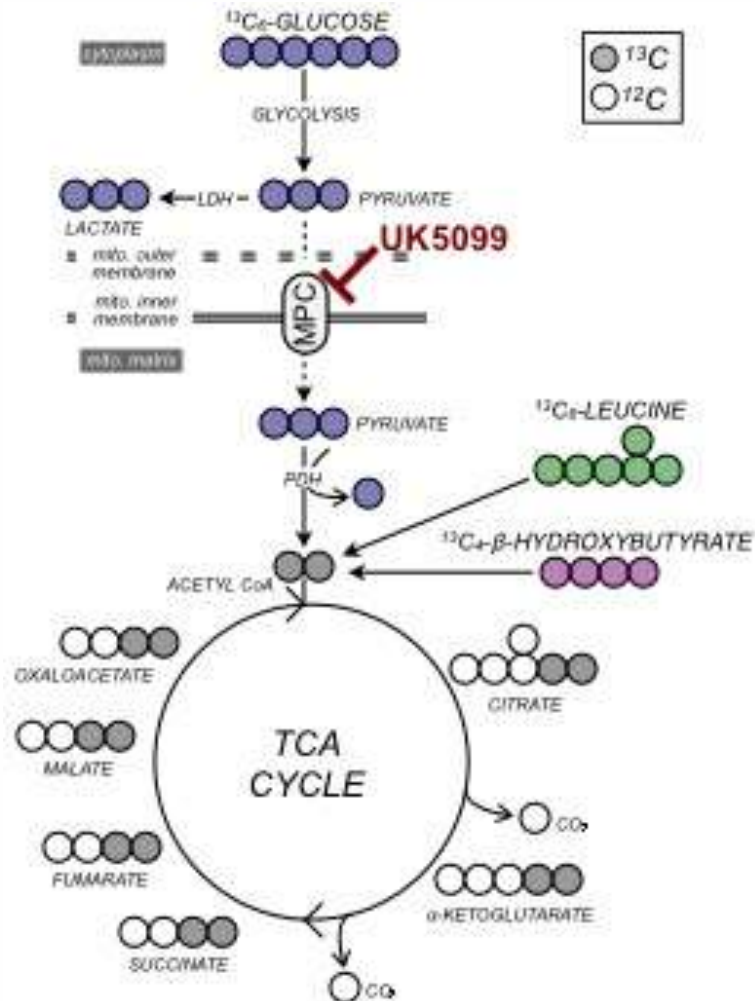
...but stable isotope tracing shows no increase in oxidation of these substrates in response to UK5099



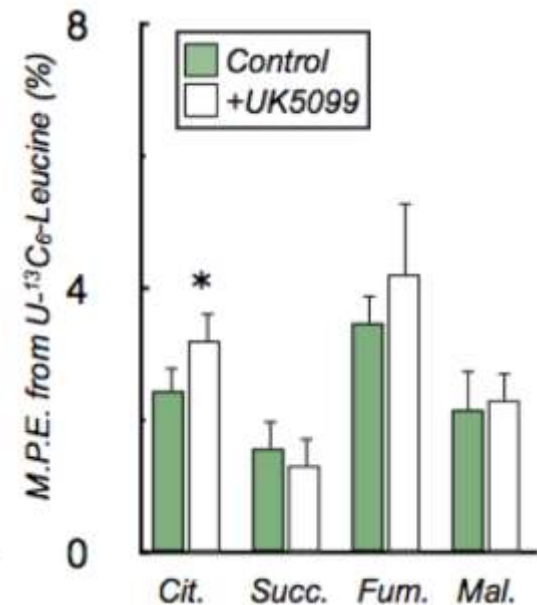
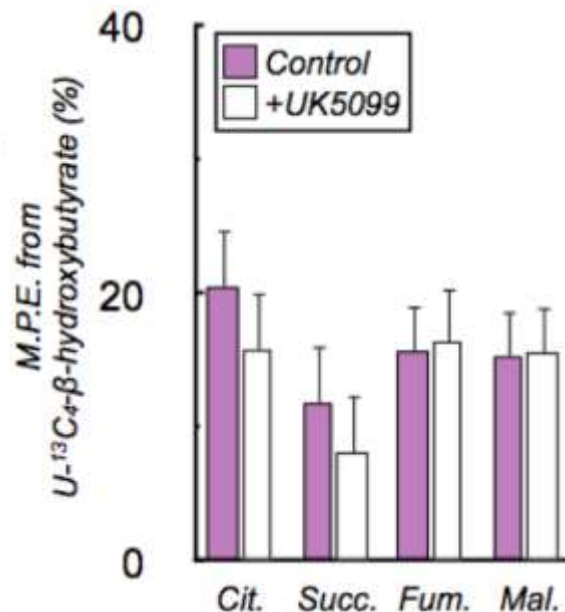
In rich culture medium we see no increase in TCA enrichment from ketone bodies (left, pink) or leucine (right, green) upon MPC inhibition over 24 hr.



...but stable isotope tracing shows no increase in oxidation of these substrates in response to UK5099

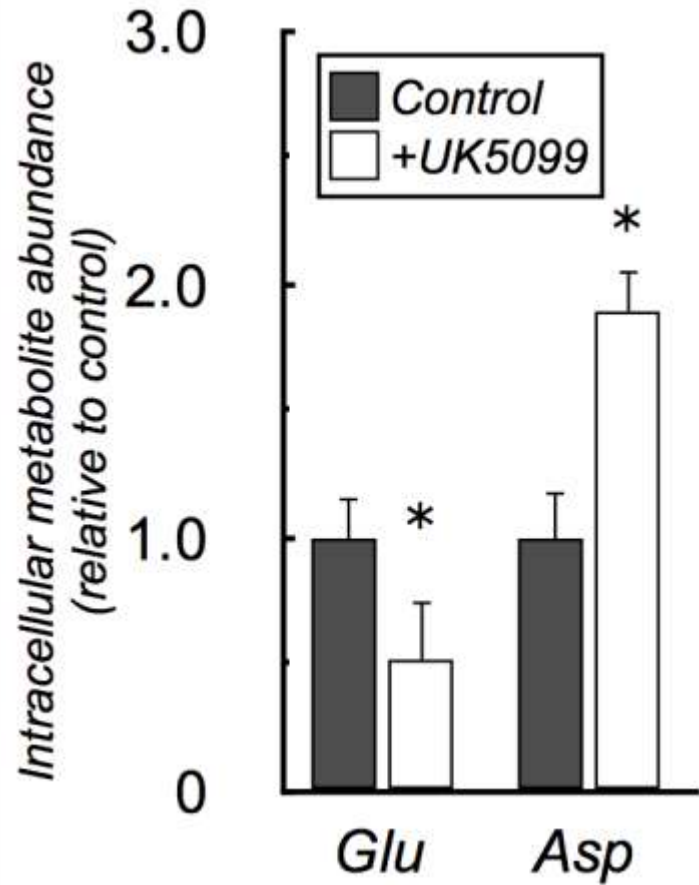
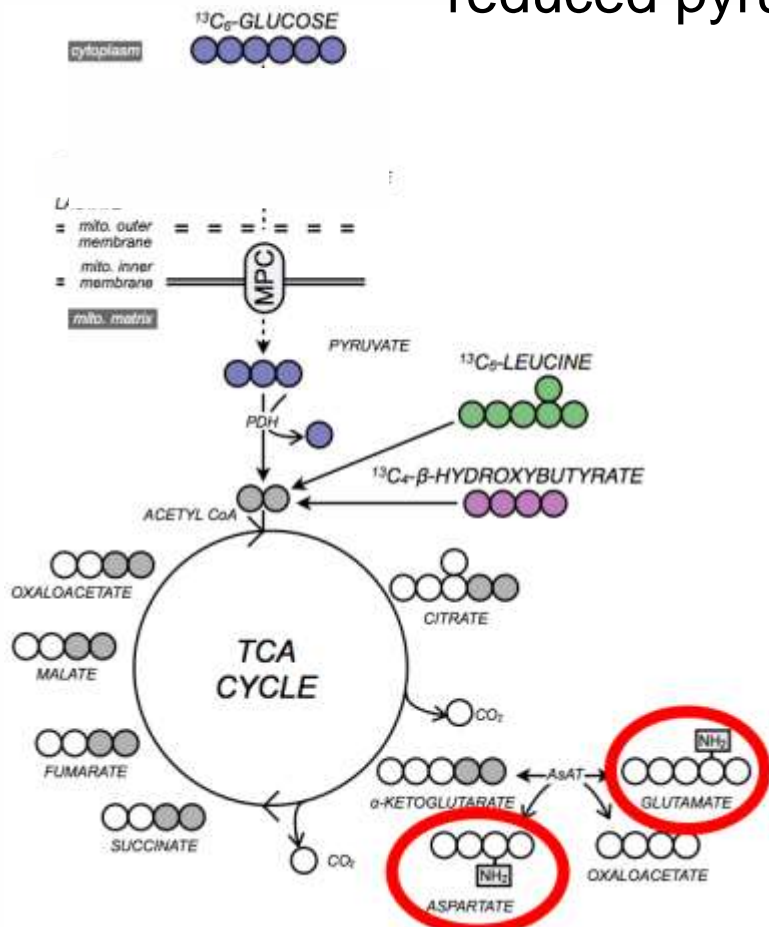


WE MISSED SOMETHING



Searching for the protective mechanism of UK5099

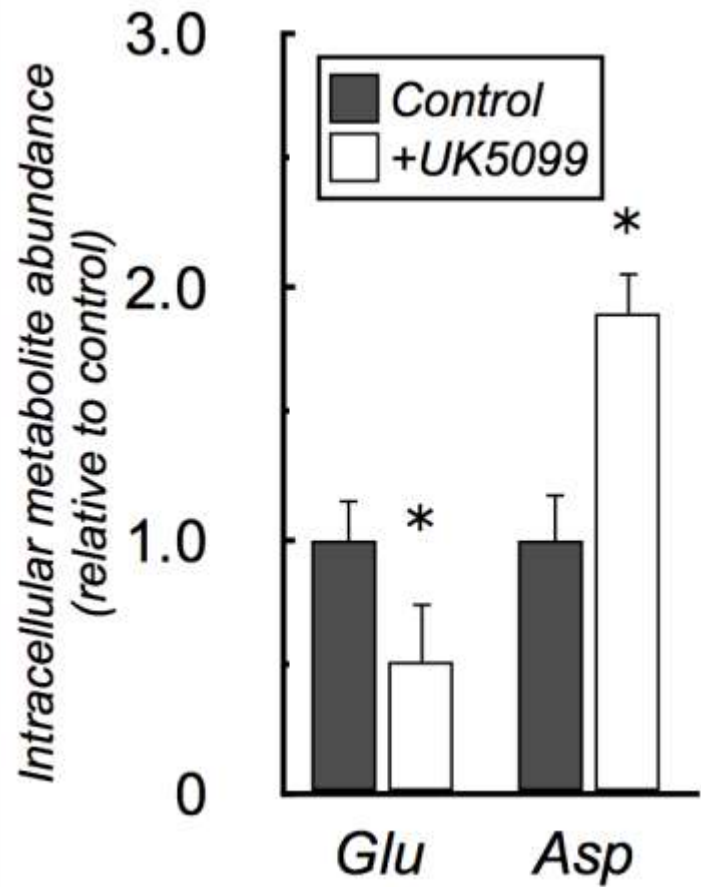
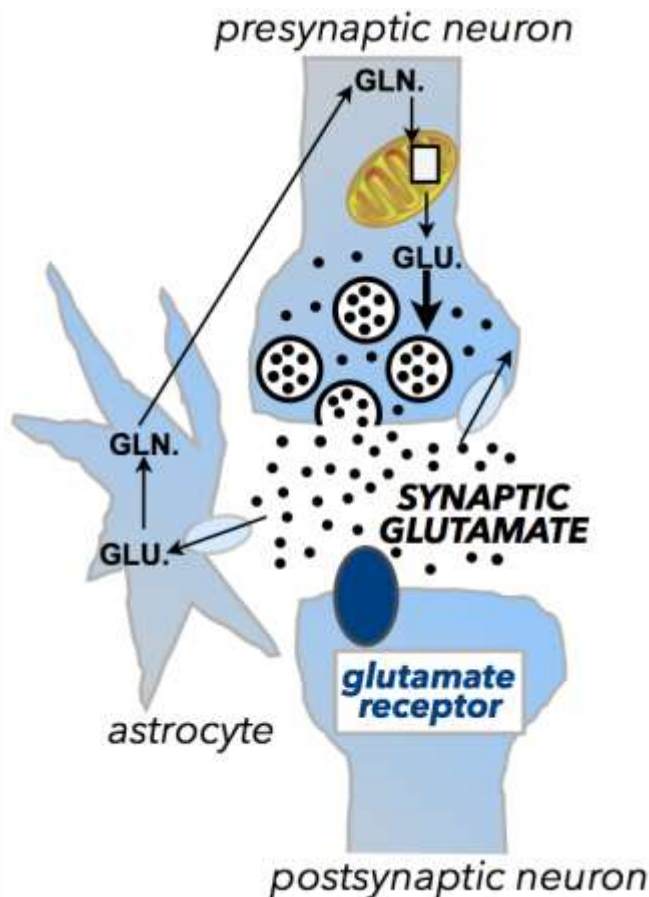
Amino acid levels suggest increased neuronal glutamate oxidation upon reduced pyruvate uptake



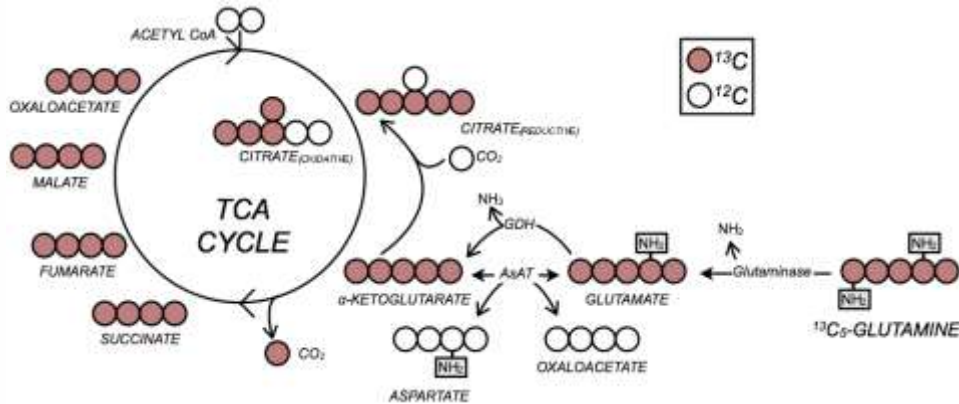
For Research Use Only. Not for use in diagnostic procedures.

Searching for the protective mechanism of UK5099

Glutamate is generally not considered a metabolic substrate due to its central (and presumably compartmentalized) role in neurotransmission

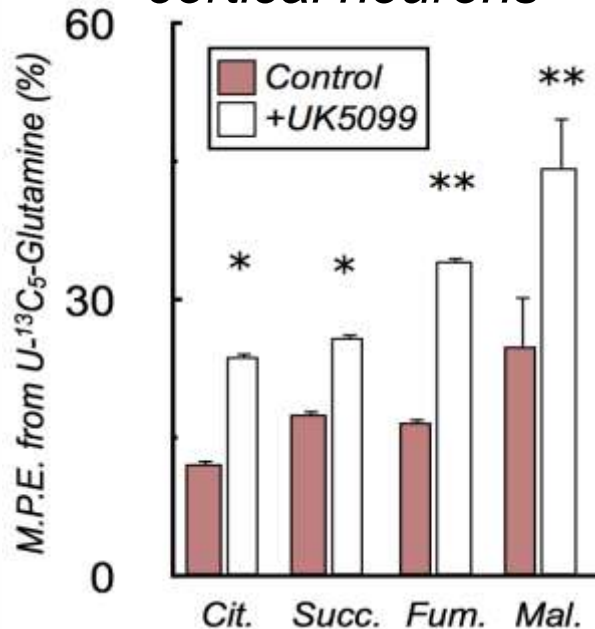


UK5099 selectively increases glutamate oxidation



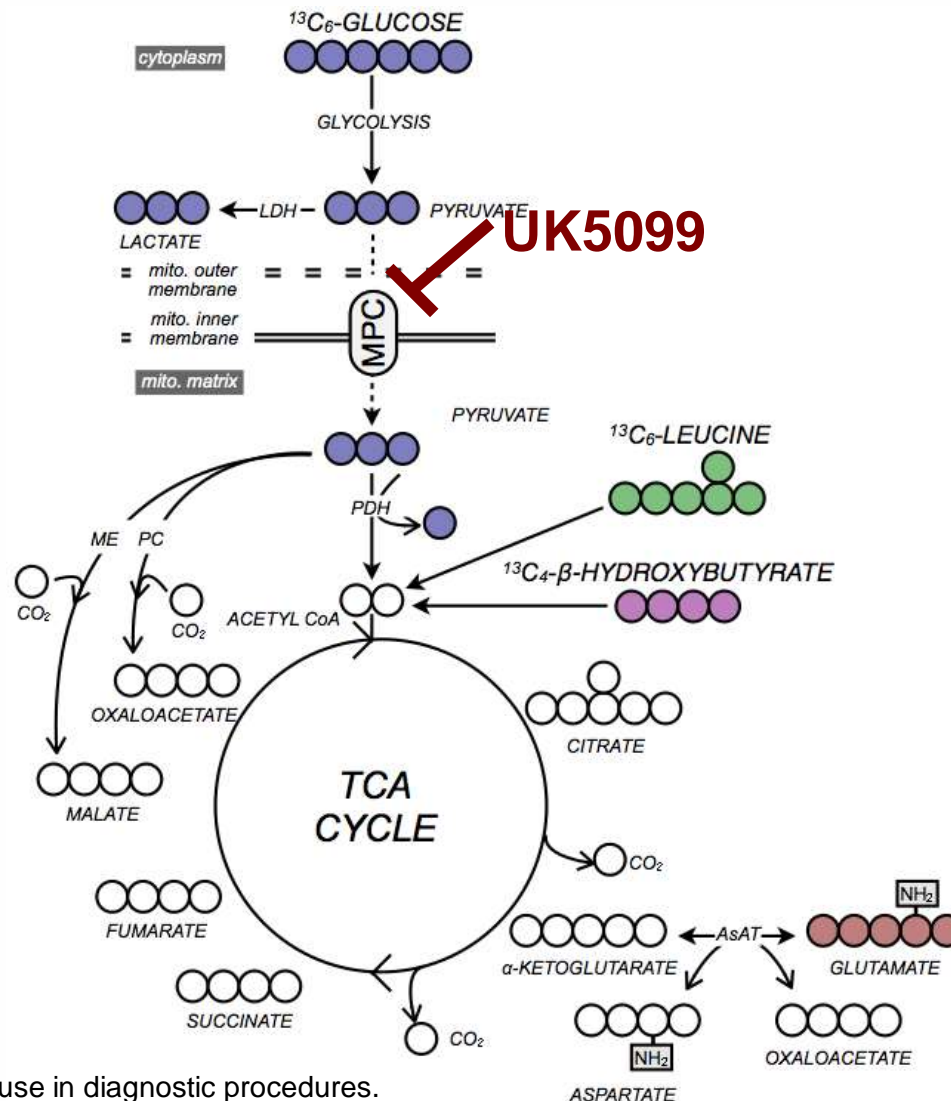
Stable isotope tracing reveals increased glutamine/glutamate oxidation in cells and organotypic slices

cortical neurons



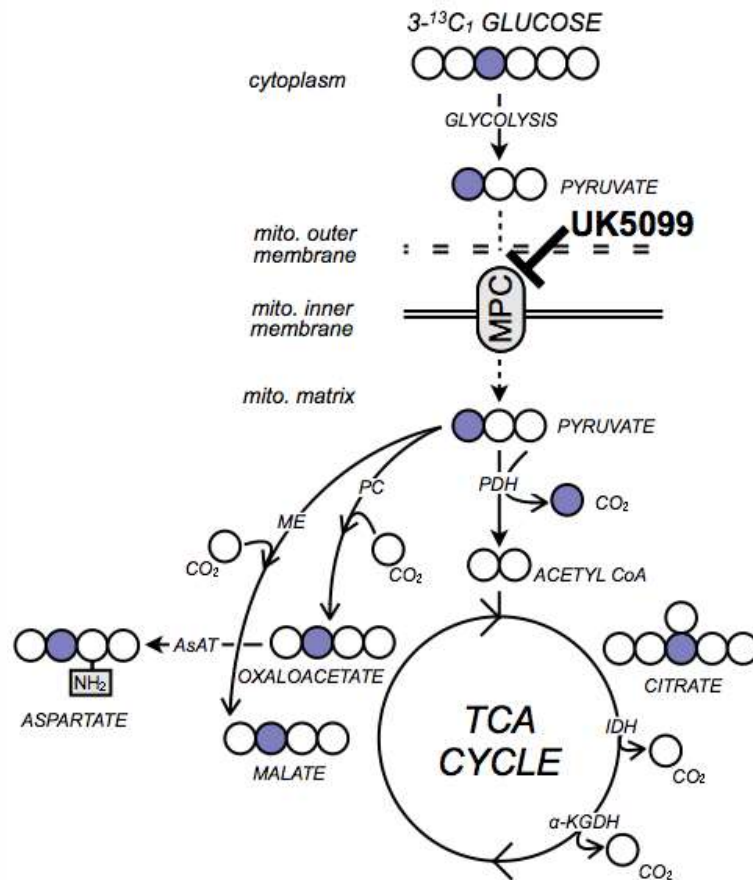
For Research Use Only. Not for use in diagnostic procedures.

Anaplerosis may explain why glutamate oxidation is selectively increased upon MPC inhibition



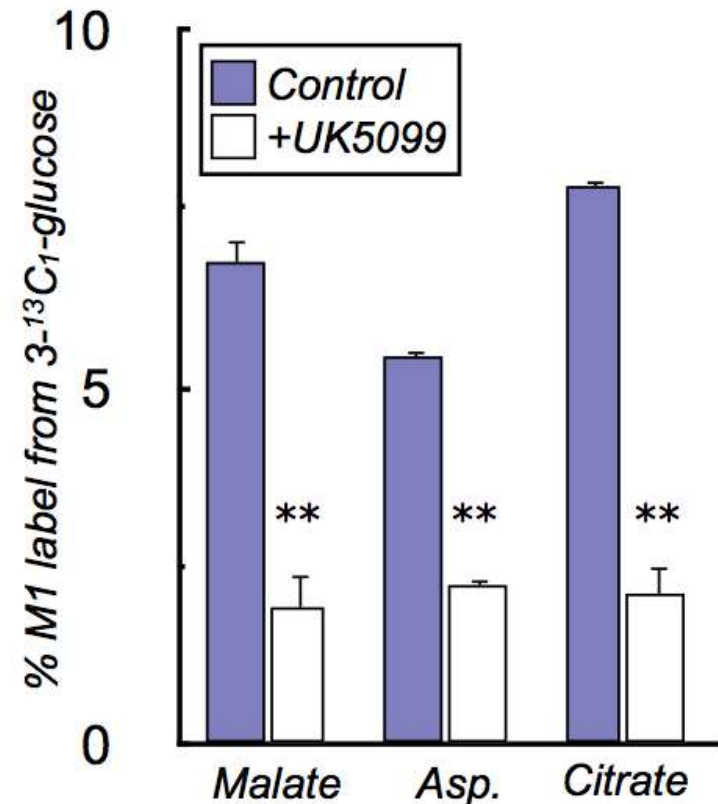
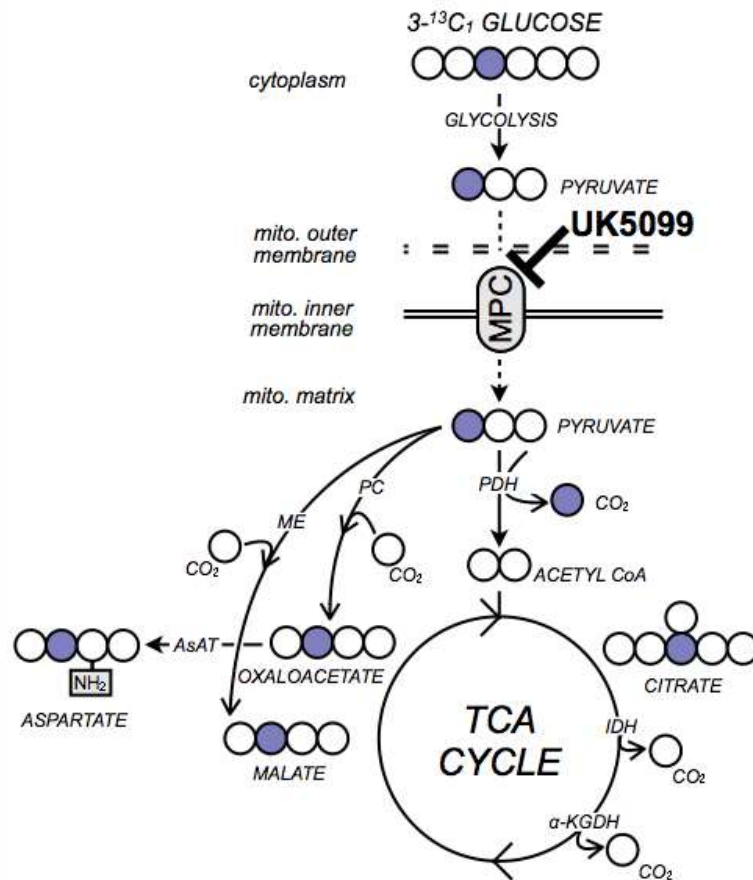
Anaplerosis may explain why glutamate oxidation is selectively increased upon MPC inhibition

A 3-¹³C₁ glucose tracer can be used to measure CO₂-fixing reactions from glucose that replenish the TCA cycle



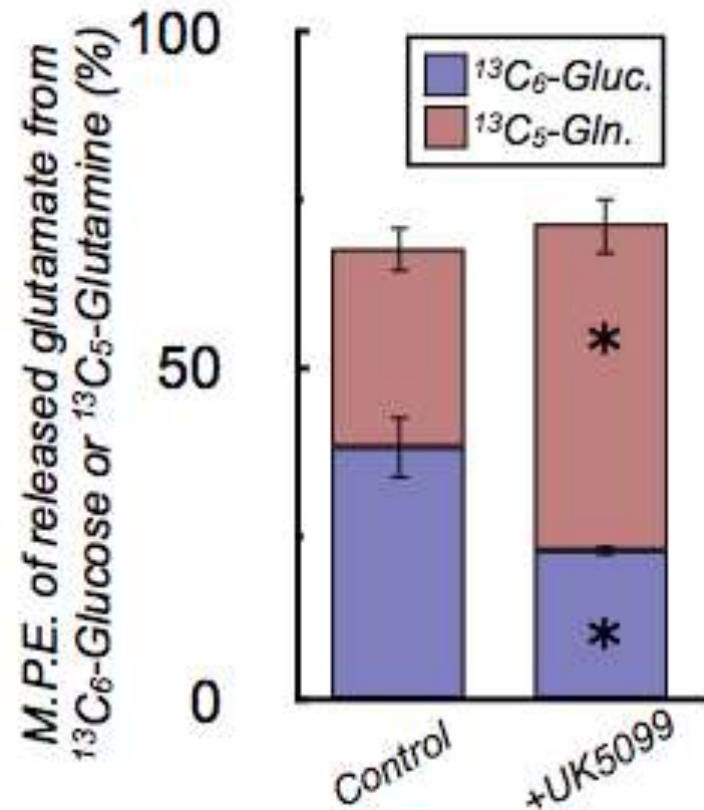
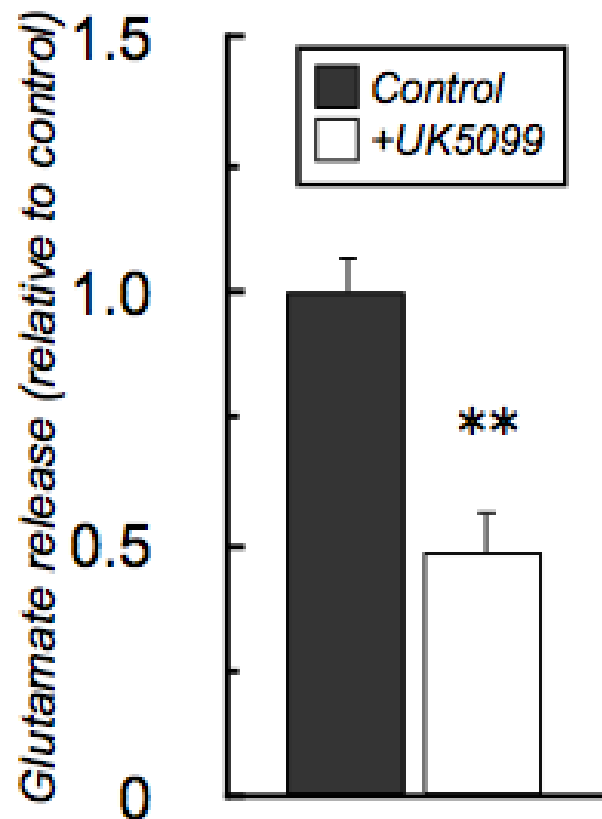
Anaplerosis may explain why glutamate oxidation is selectively increased upon MPC inhibition

A 3- $^{13}\text{C}_1$ glucose tracer can be used to measure CO_2 -fixing reactions from glucose that replenish the TCA cycle



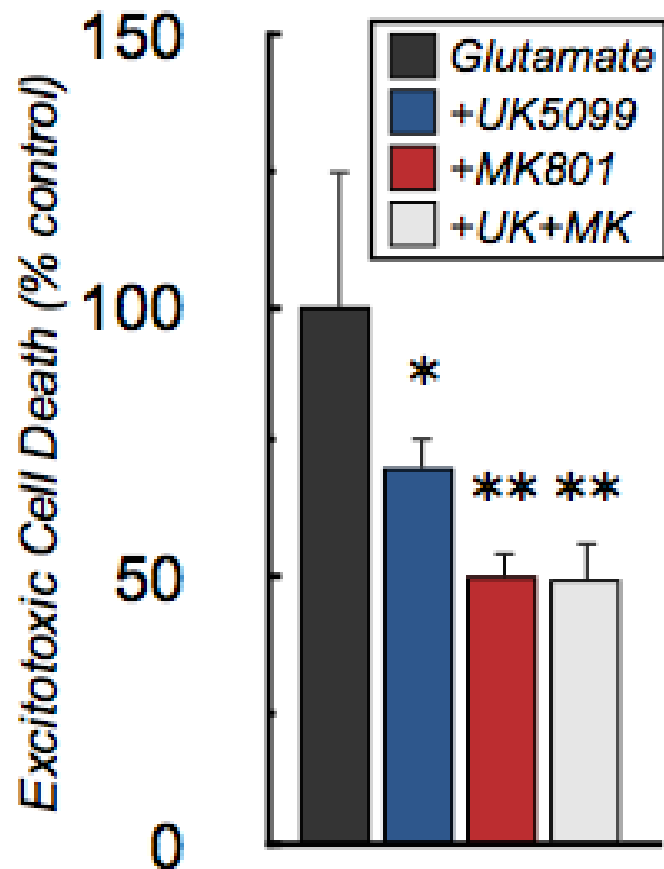
The glutamate mobilized for oxidation is not discrete from the pool apportioned for evoked release

Glutamate released by depolarization from veratridine (prevents Na^+ -channel closure) and ouabain (Na^+/K^+ -ATPase inhibitor)



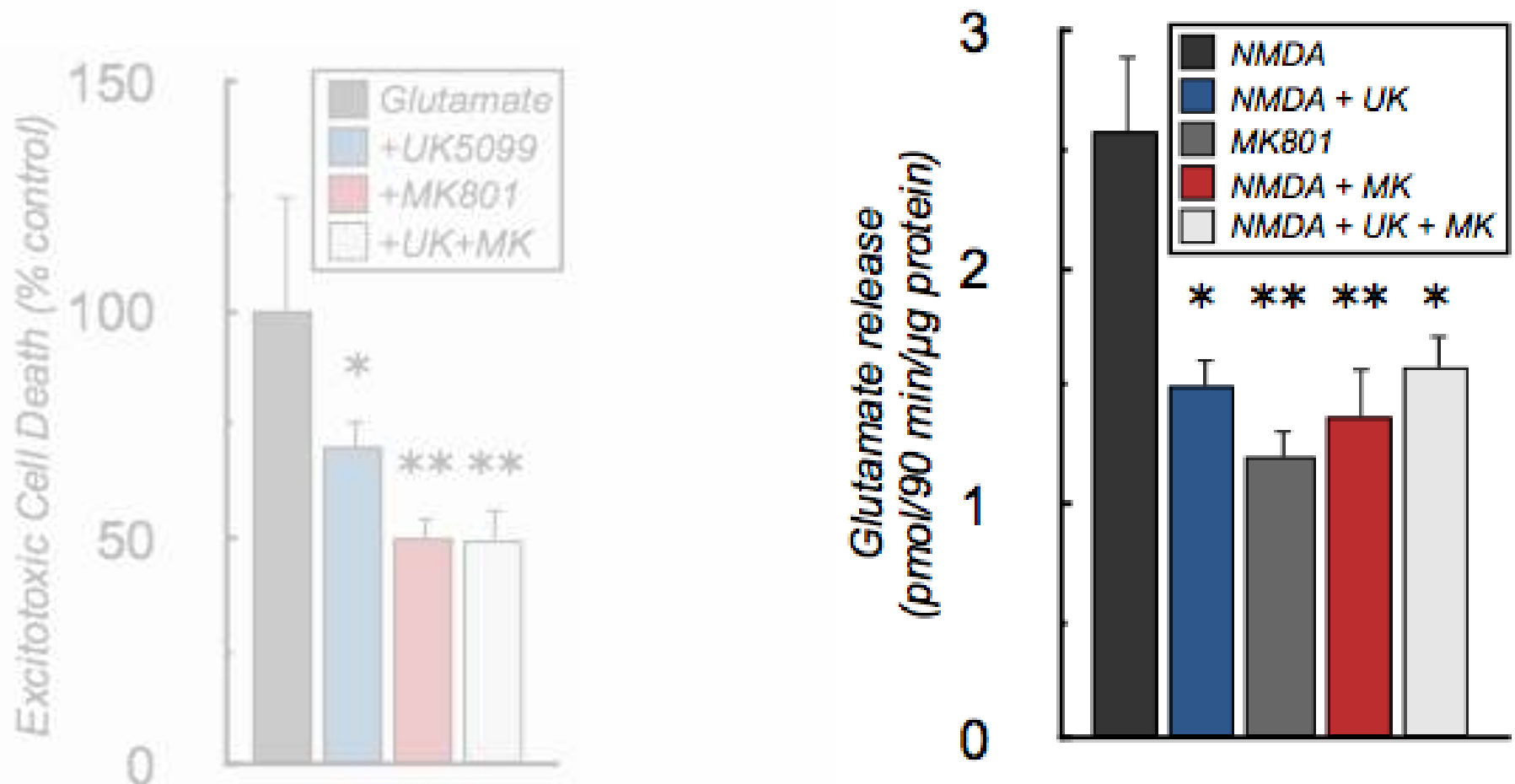
Can attenuated vesicular release explain the protection from excitotoxic injury?

Limiting depolarization-induced glutamate release could minimize the positive-feedback, excitotoxic cascade



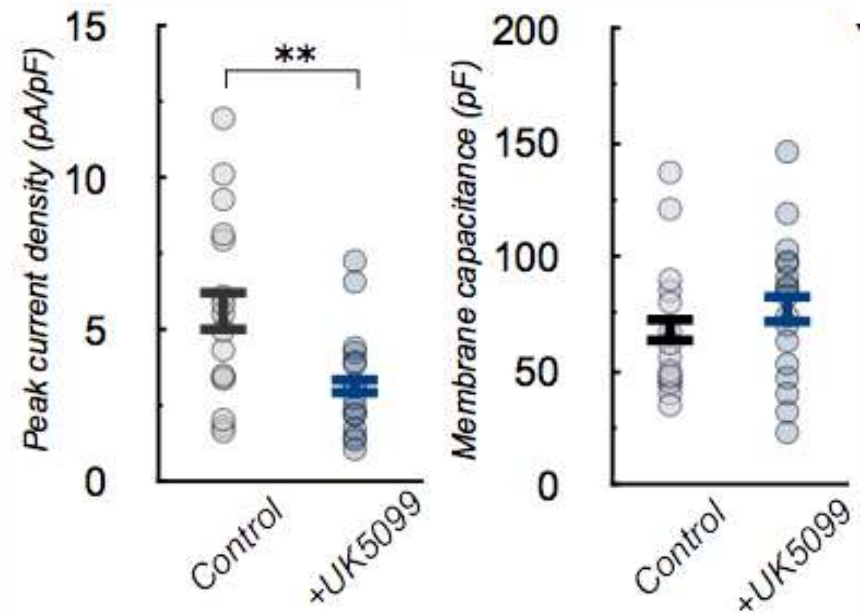
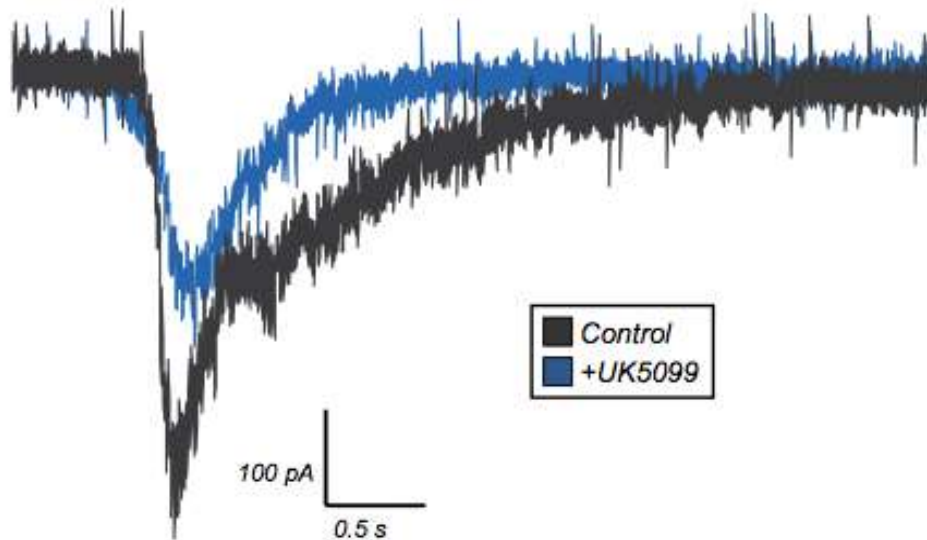
Can attenuated vesicular release explain the protection from excitotoxic injury?

Depolarization induced with a concentration & duration of NMDA that does not compromise viability

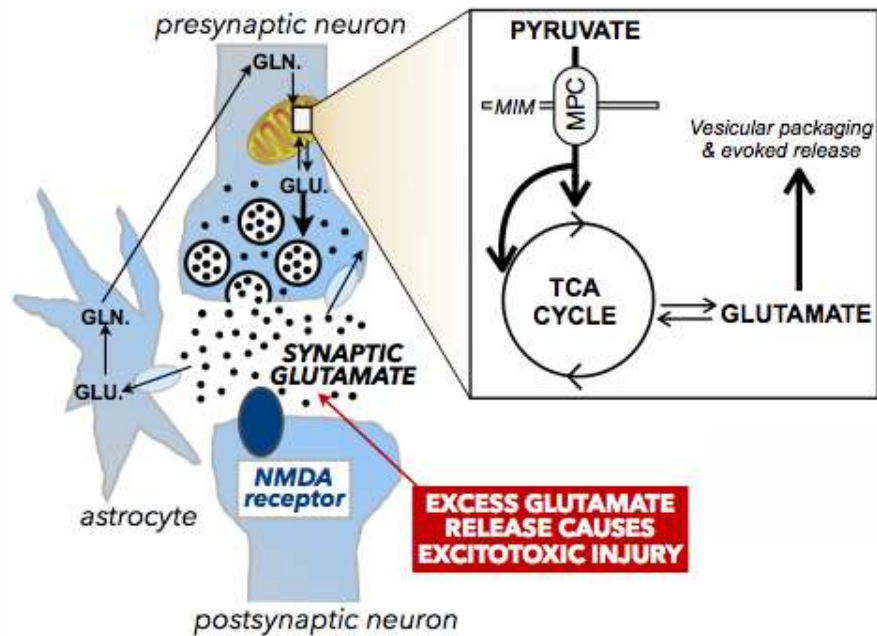


Pyruvate carrier inhibition lowers synaptic glutamate release

UK5099 attenuates excitatory post-synaptic currents in primary mouse pyramidal neurons



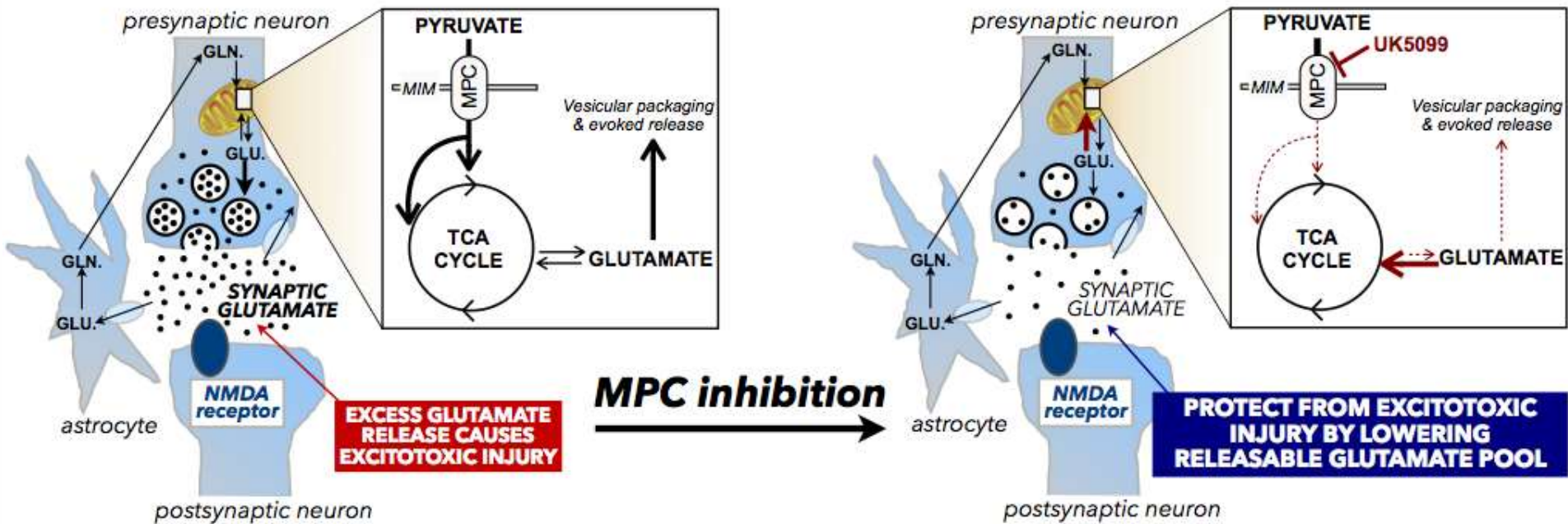
Our model



Divakaruni...and Murphy (2017) J. Cell Biol.

For Research Use Only. Not for use in diagnostic procedures.

Our model



Divakaruni...and Murphy (2017) J. Cell Biol.

For Research Use Only. Not for use in diagnostic procedures.

Acknowledgements



Anne Murphy

Sandy Wiley Brenda Bloodgood

Alex Andreyev Kelly Martyniuk



Lynn Raymond
Caodu Buren

FUNDING

NIH (NS087611, DK05441, DK81298)

AGILENT TECHNOLOGIES

TEVA PHARMACEUTICALS



Christian Metallo

Martina Wallace

Nate Vacanti



Agilent Technologies