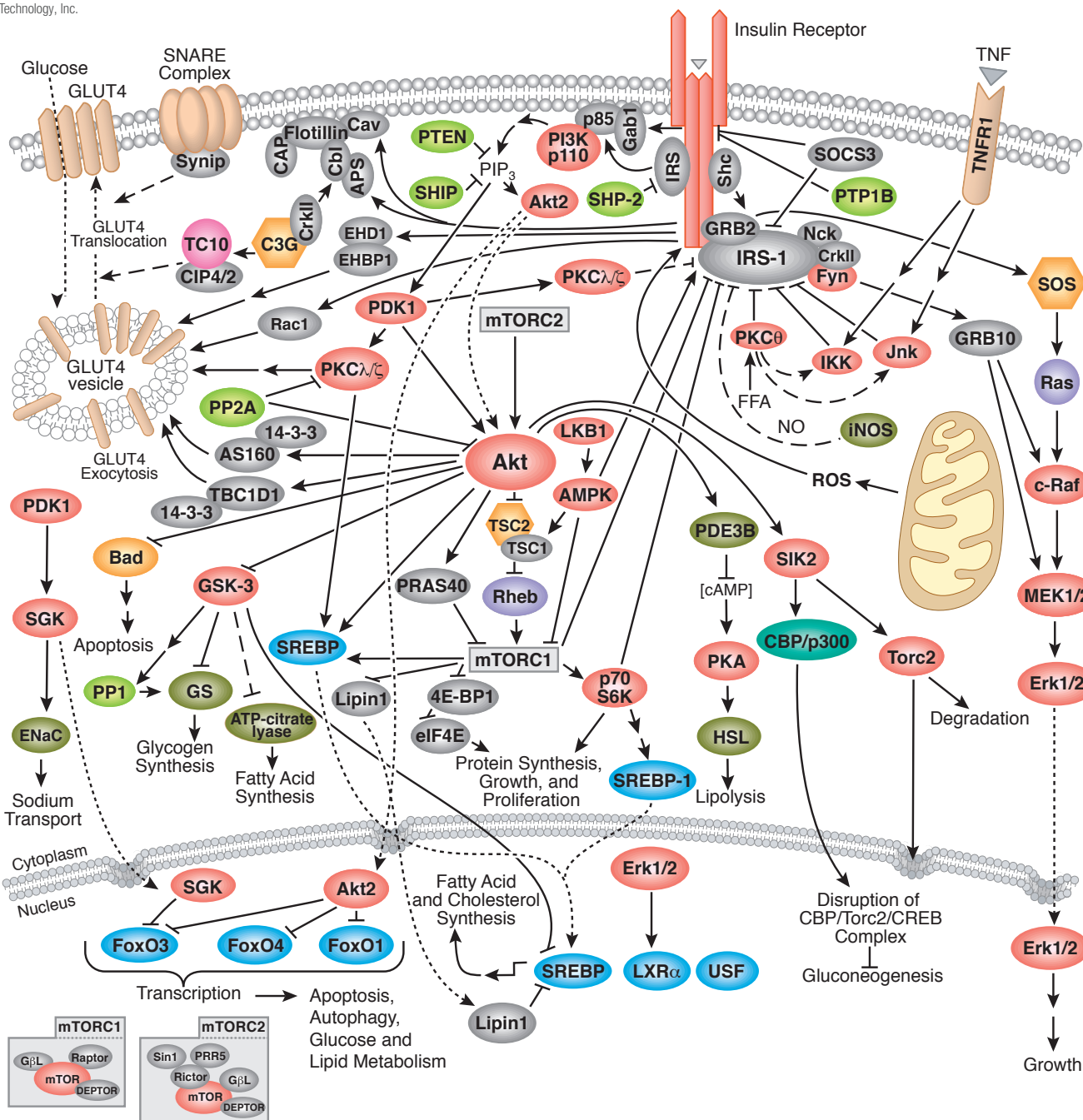


Cellular Metabolism

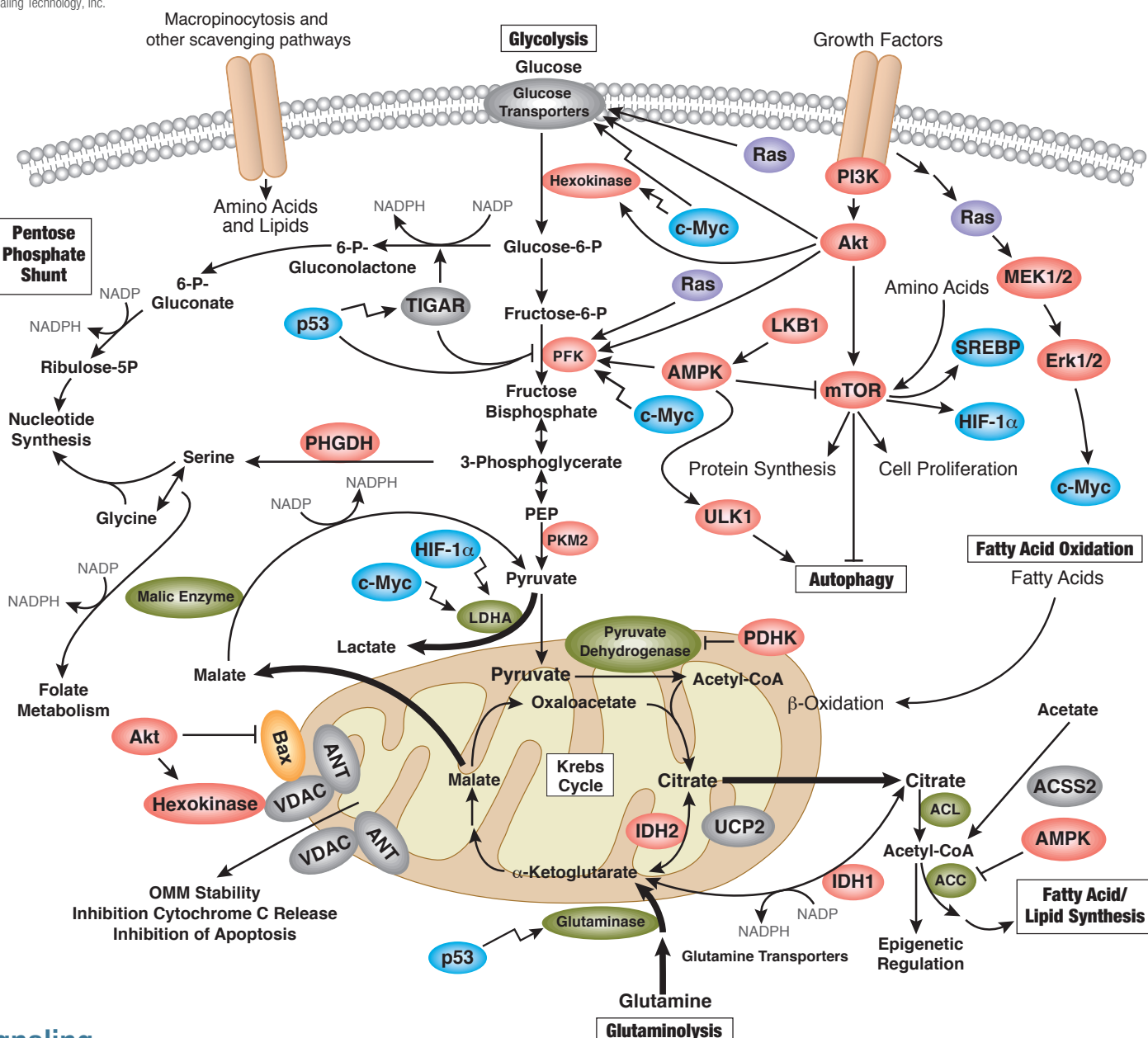
Insulin Receptor Signaling

© 2003–2015 Cell Signaling Technology, Inc.



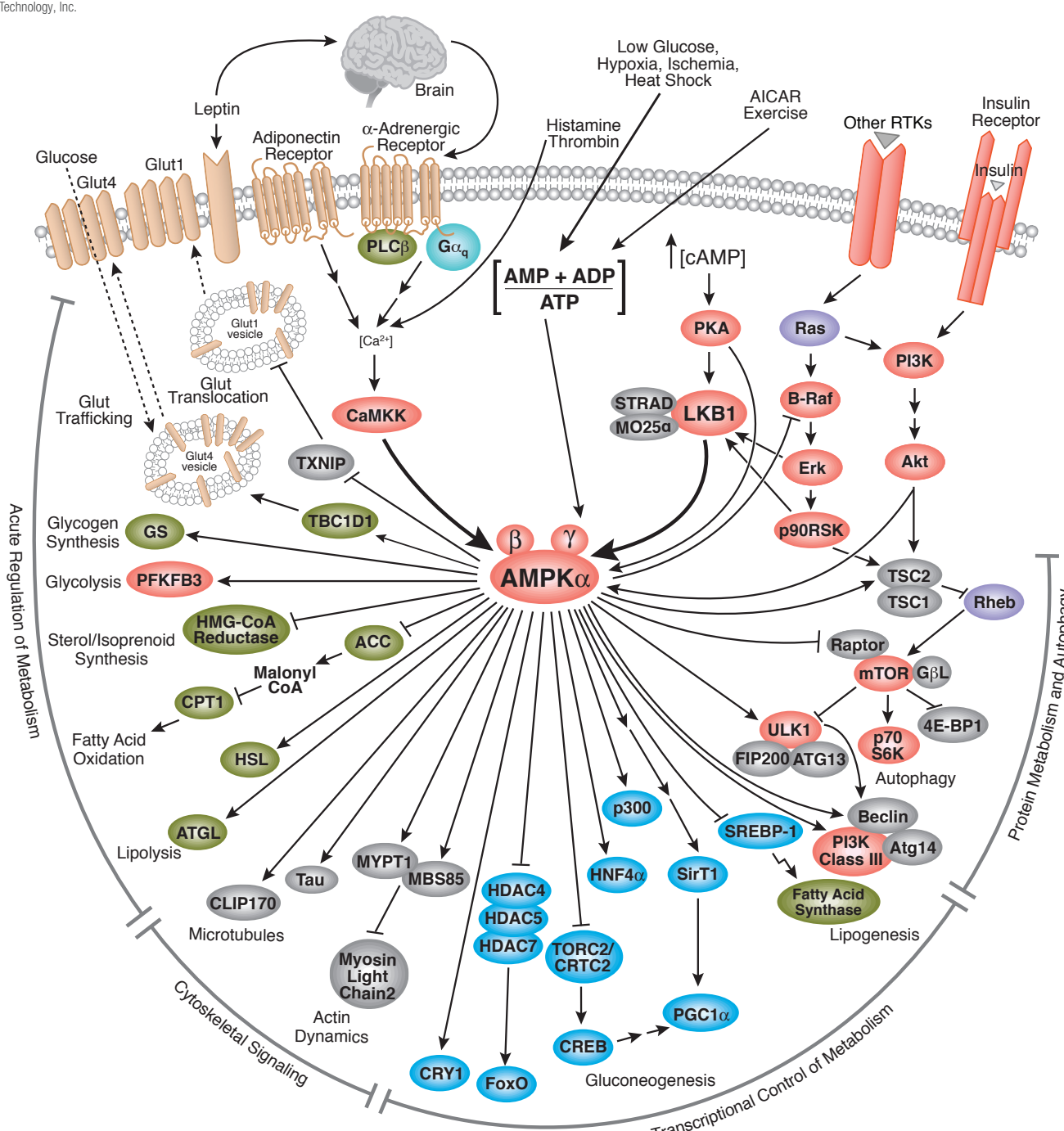
Warburg Effect

© 2010–2015 Cell Signaling Technology, Inc.



AMPK Signaling

© 2008–2015 Cell Signaling Technology, Inc.



→ Direct Stimulatory Modification	→ Multistep Inhibitory Modification	→ Separation of Subunits or Cleavage Products	→ Transcriptional Inhibitory Modification	○ Receptor	○ GTPase
→ Direct Inhibitory Modification	→ Tentative Stimulatory Modification	→ Joining of Subunits	○ Kinase	○ Enzyme	○ G-protein
→ Multistep Stimulatory Modification	→ Tentative Inhibitory Modification	→ Translocation	○ Phosphatase	○ pro-apoptotic	○ Acetylase
		→ Transcriptional Stimulatory Modification	○ Transcription Factor	○ pro-survival	○ Deacetylase
			○ Caspase	○ GAP/GEF	○ Ribosomal subunit

Printed in the USA on recycled paper (25% post-consumer waste fiber) using vegetable inks and processed chlorine free. © 2015 Cell Signaling Technology, Inc. Cell Signaling Technology® and CST™ are trademarks of Cell Signaling Technology, Inc.