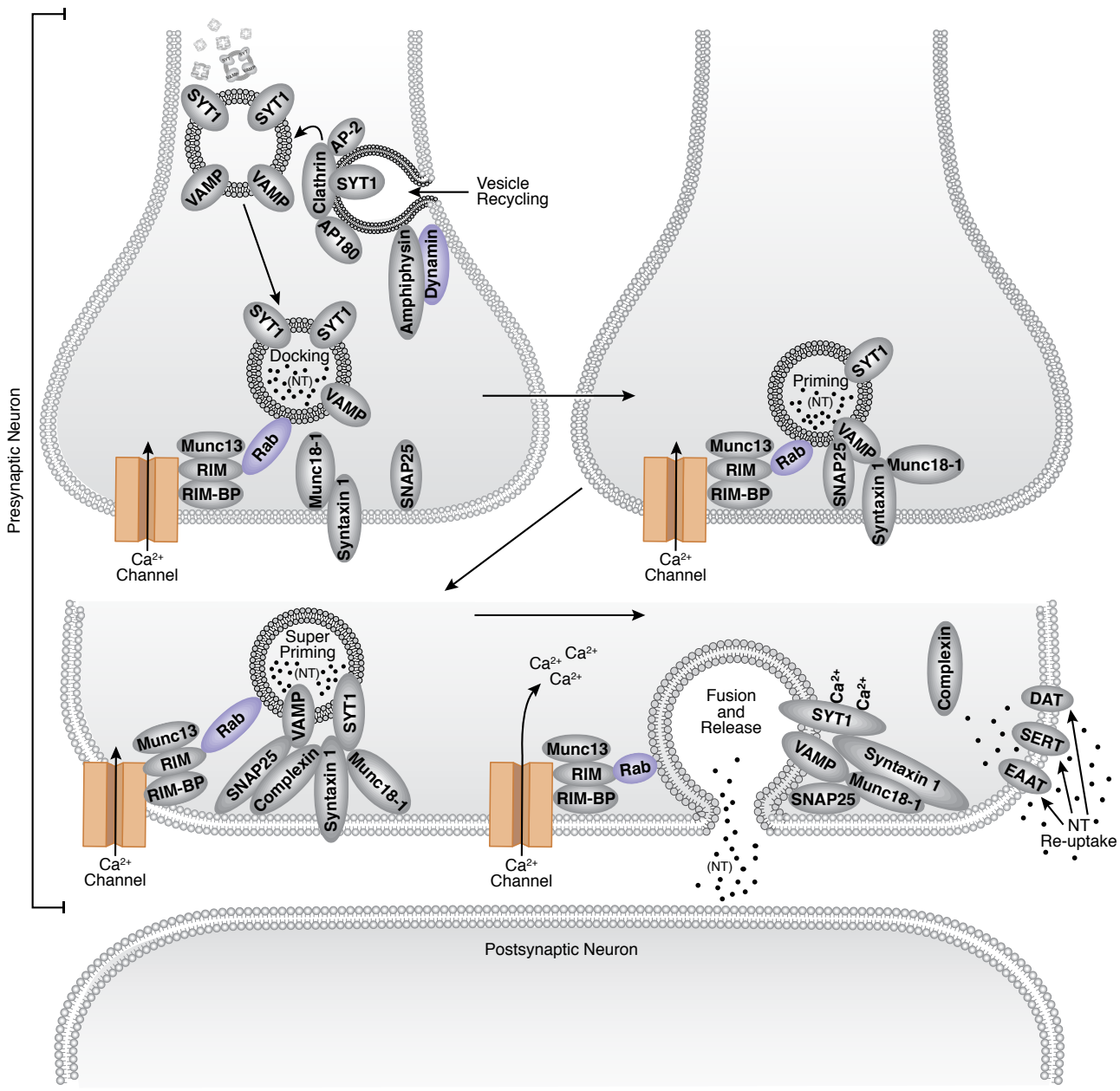


# Neuroscience

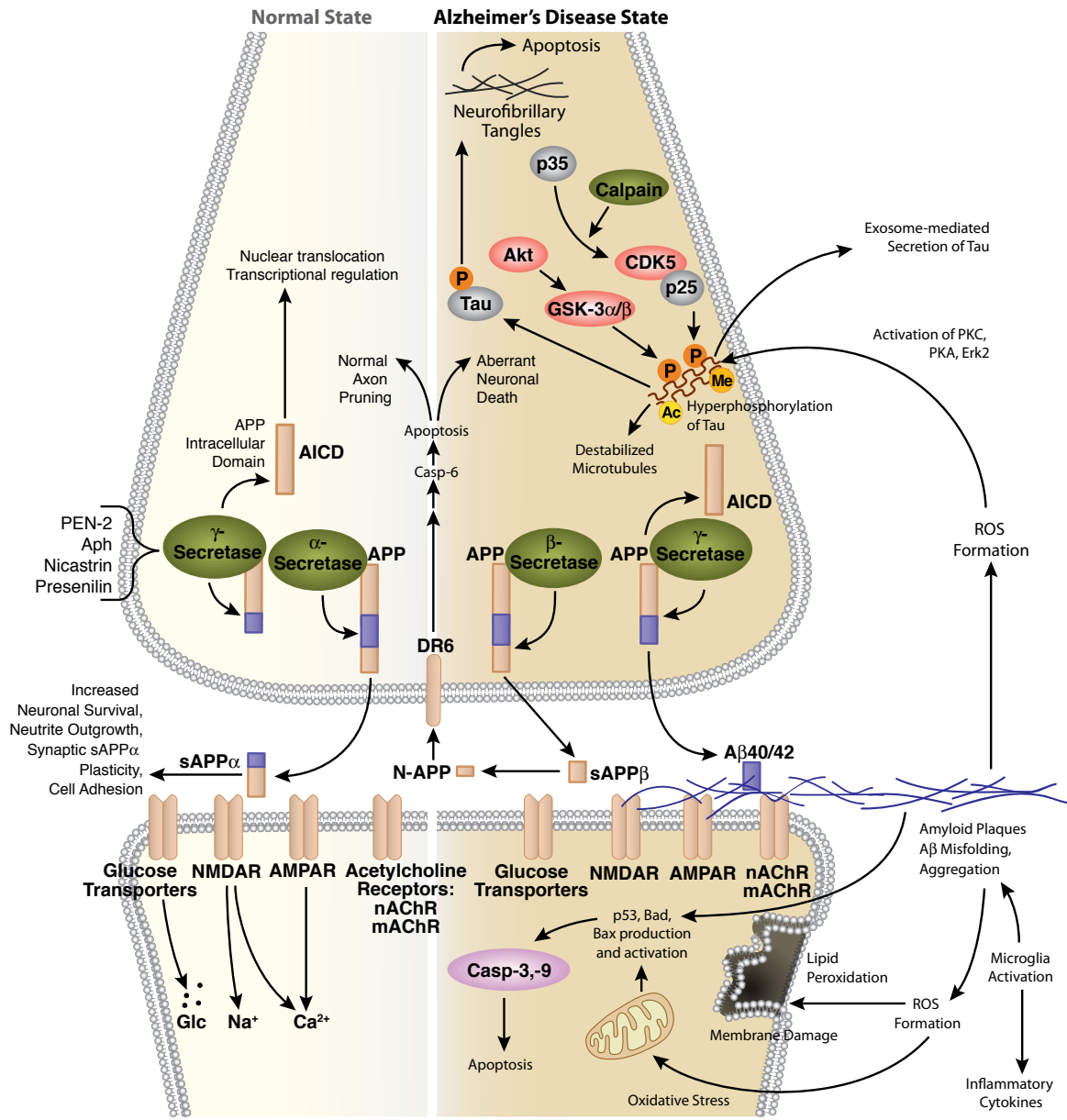
## Vesicle Trafficking in Presynaptic Neurons: Synchronous Release

© 2014–2015 Cell Signaling Technology, Inc.



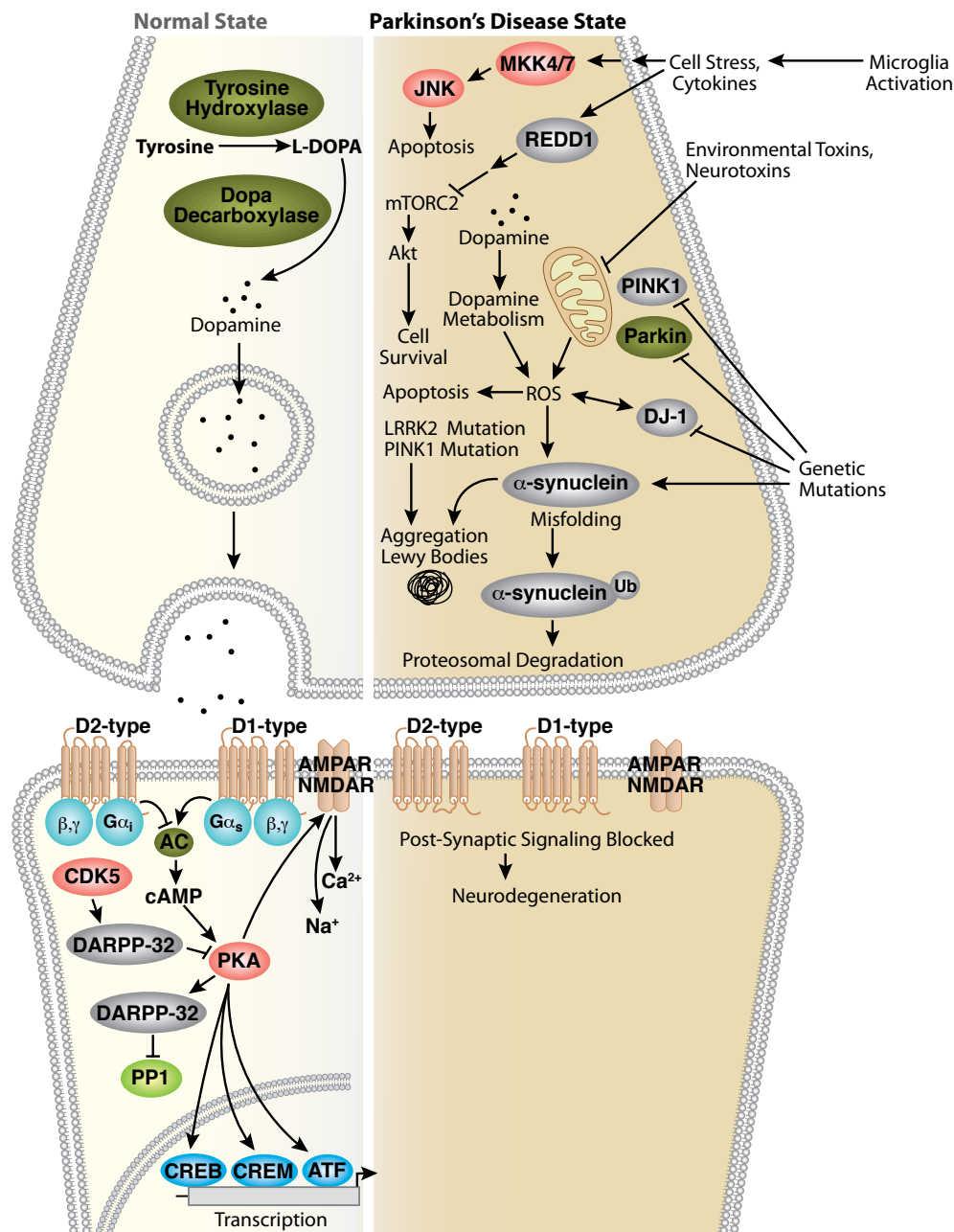
## Amyloid Plaque and Neurofibrillary Tangle Formation in Alzheimer's Disease

© 2009–2015 Cell Signaling Technology, Inc.



## Dopamine Signaling in Parkinson's Disease

© 2009–2015 Cell Signaling Technology, Inc.



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