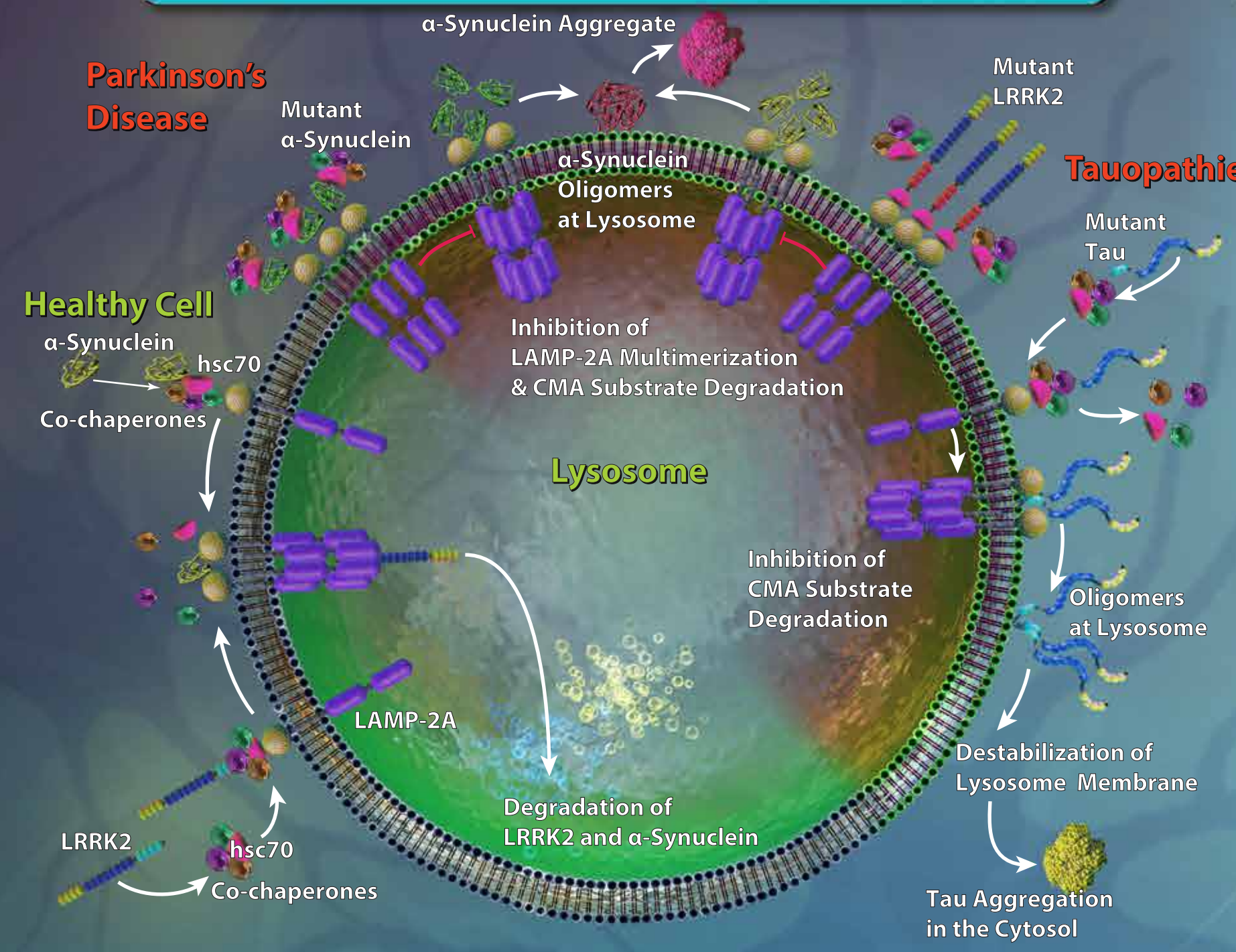
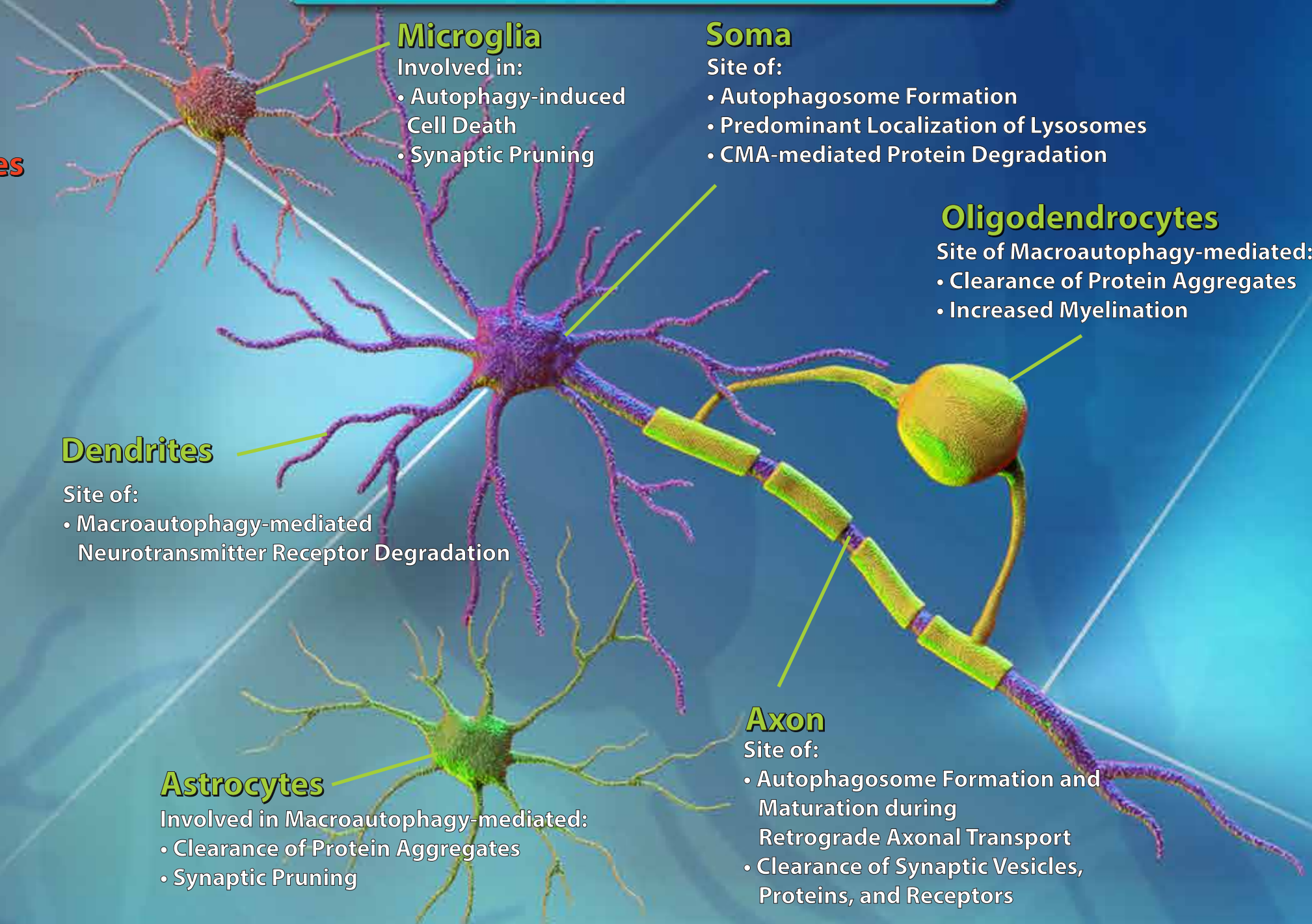


Autophagy

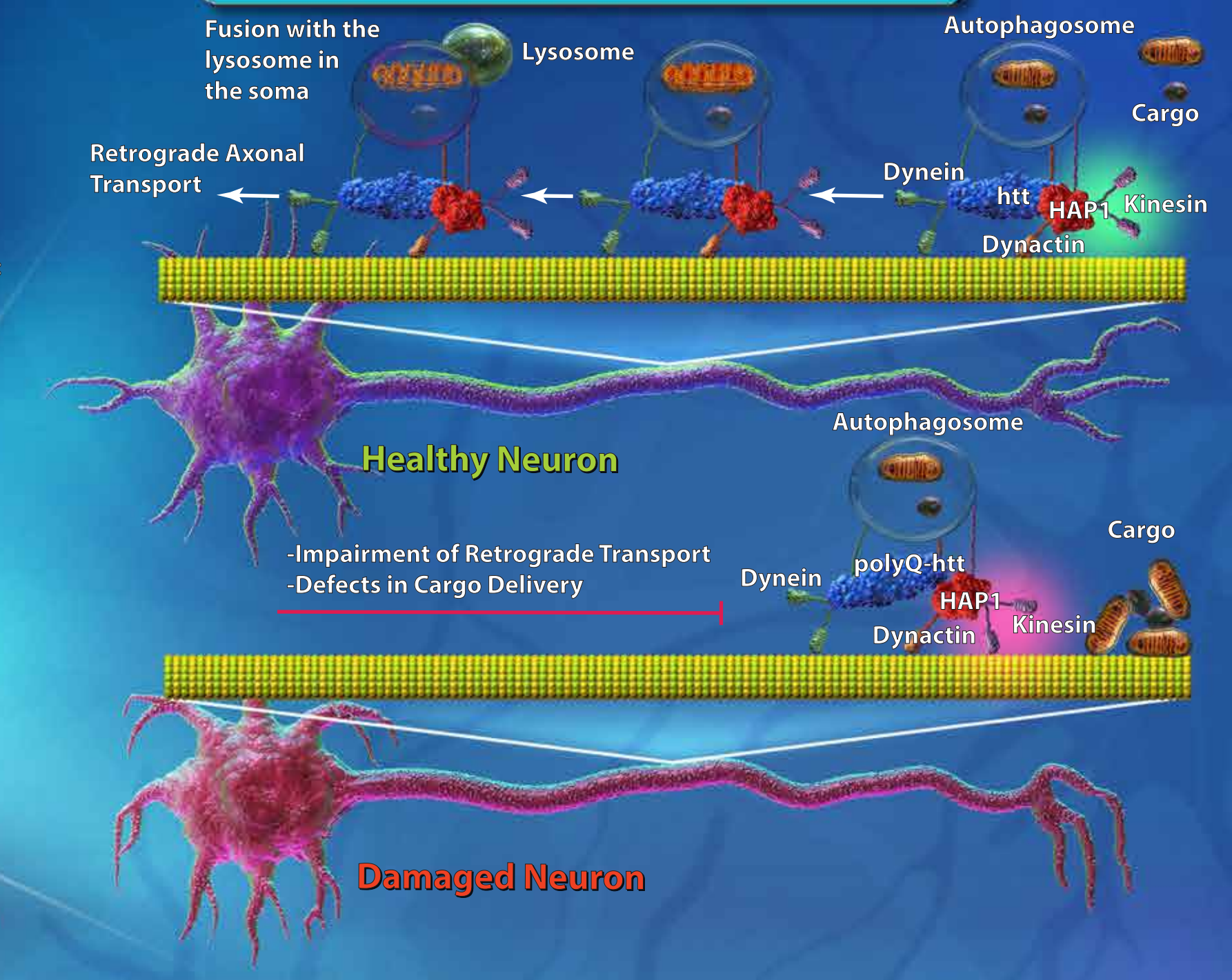
Chaperone-mediated Autophagy (ATG Independent)



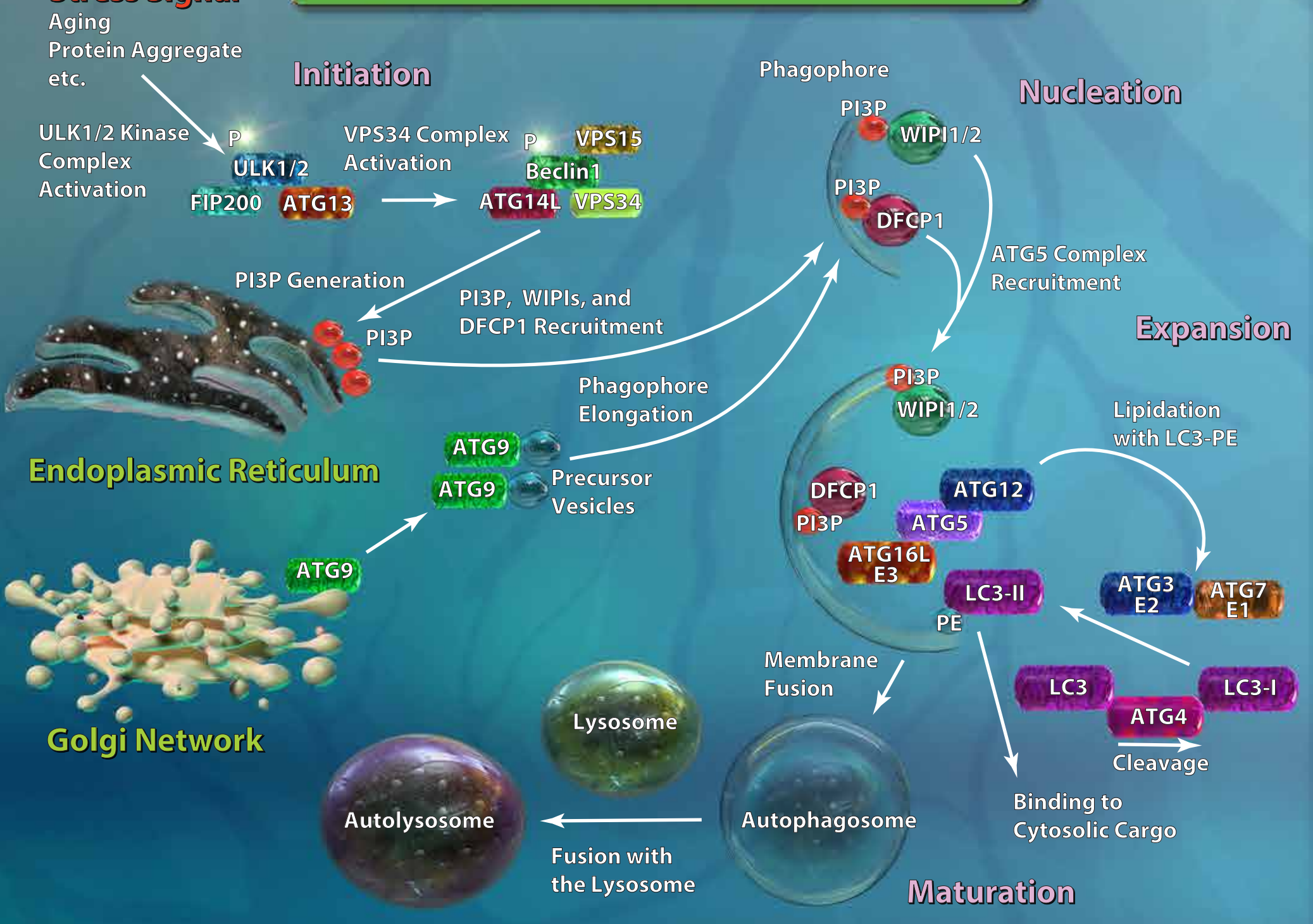
Autophagy Events in the Brain



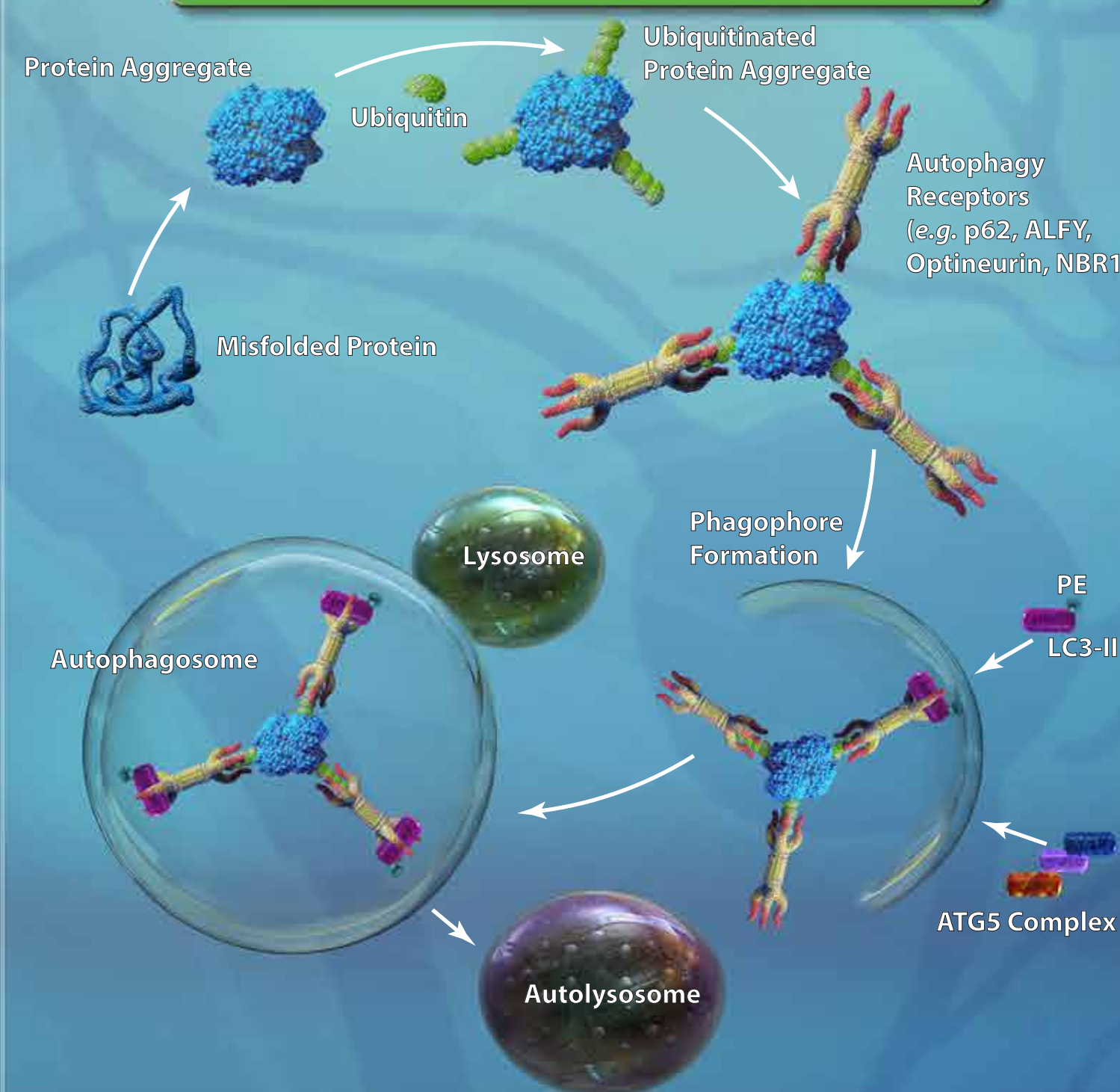
Autophagy in Axonal Transport



Macroautophagy Process (ATG Dependent)



Aggrephagy (Selective Macroautophagy)



Mitophagy (Selective Macroautophagy)

