

# Protein Recycling, TPD, and Neurodegenerative Disease



Cells have two main strategies for degrading and recycling proteins:

- 1) The ubiquitin (Ub) proteasome system (UPS)
- 2) Autophagy.

However, faults in proteins or these degradation systems can lead to neurodegenerative diseases, such as Parkinson's, Alzheimer's, and Huntington's disease.

Targeted protein degradation (TPD) is a therapeutic strategy that hijacks protein degradation pathways, which could potentially treat neurodegenerative diseases.

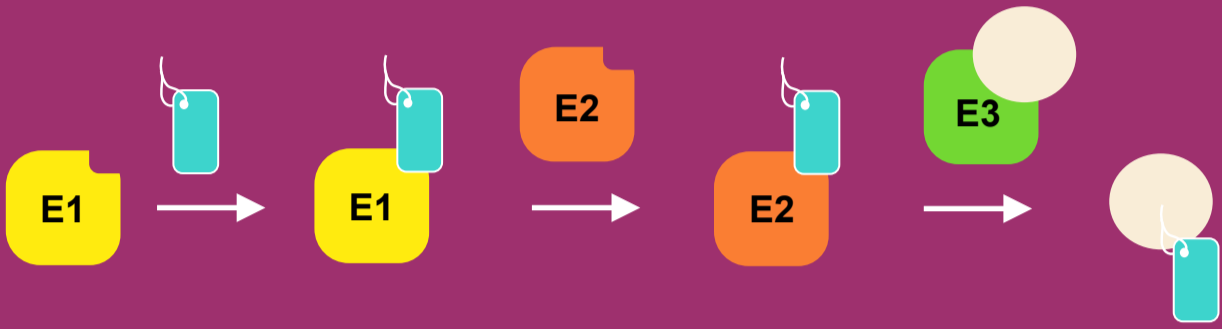


## Protein Recycling Centre

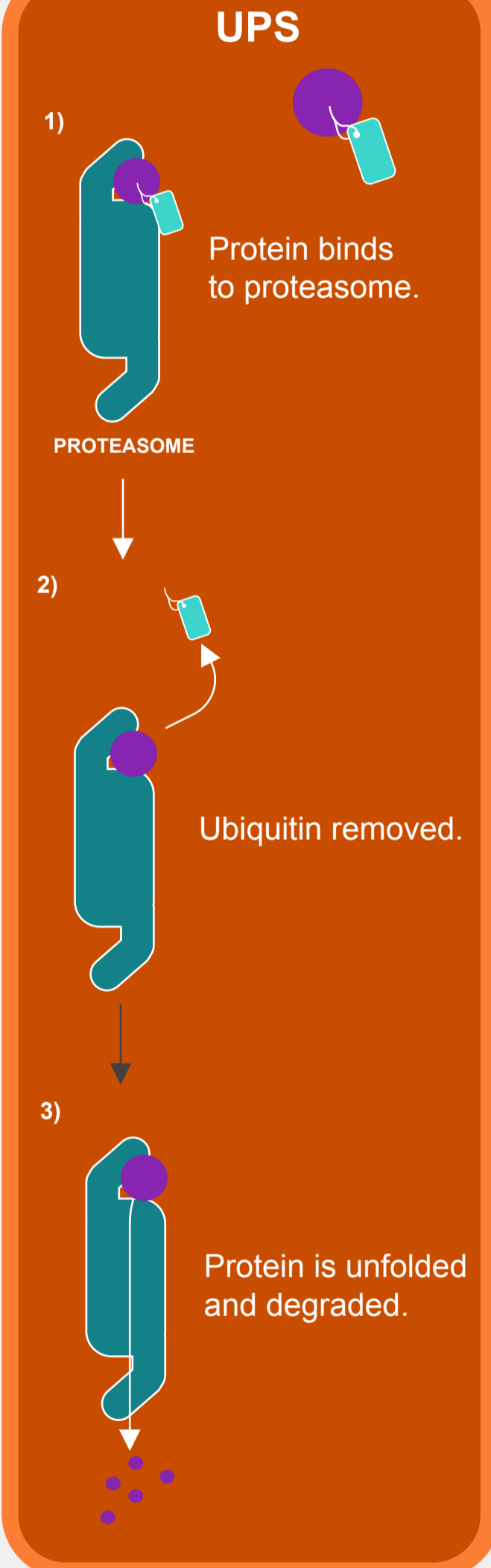
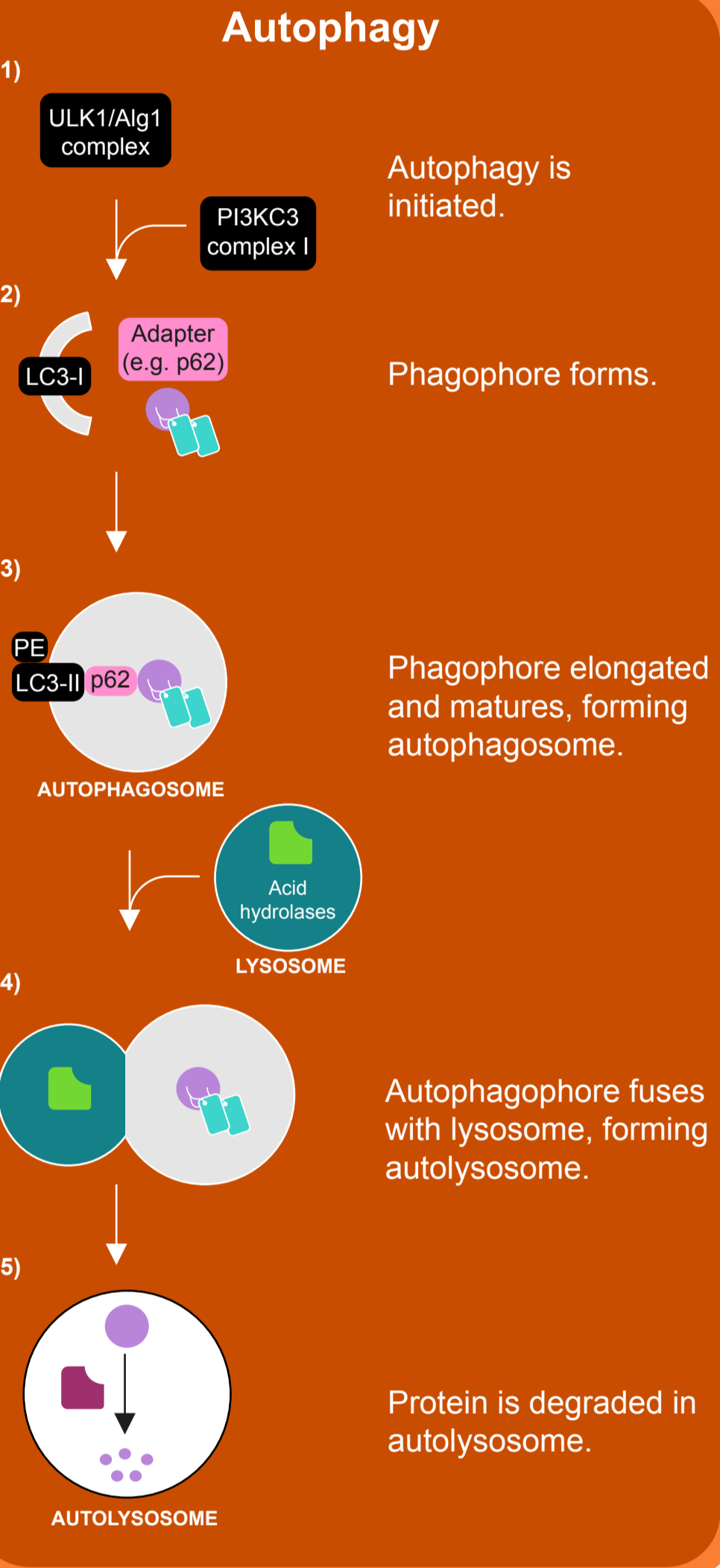


**Sorting Station:**  
Ubiquitination of Proteins

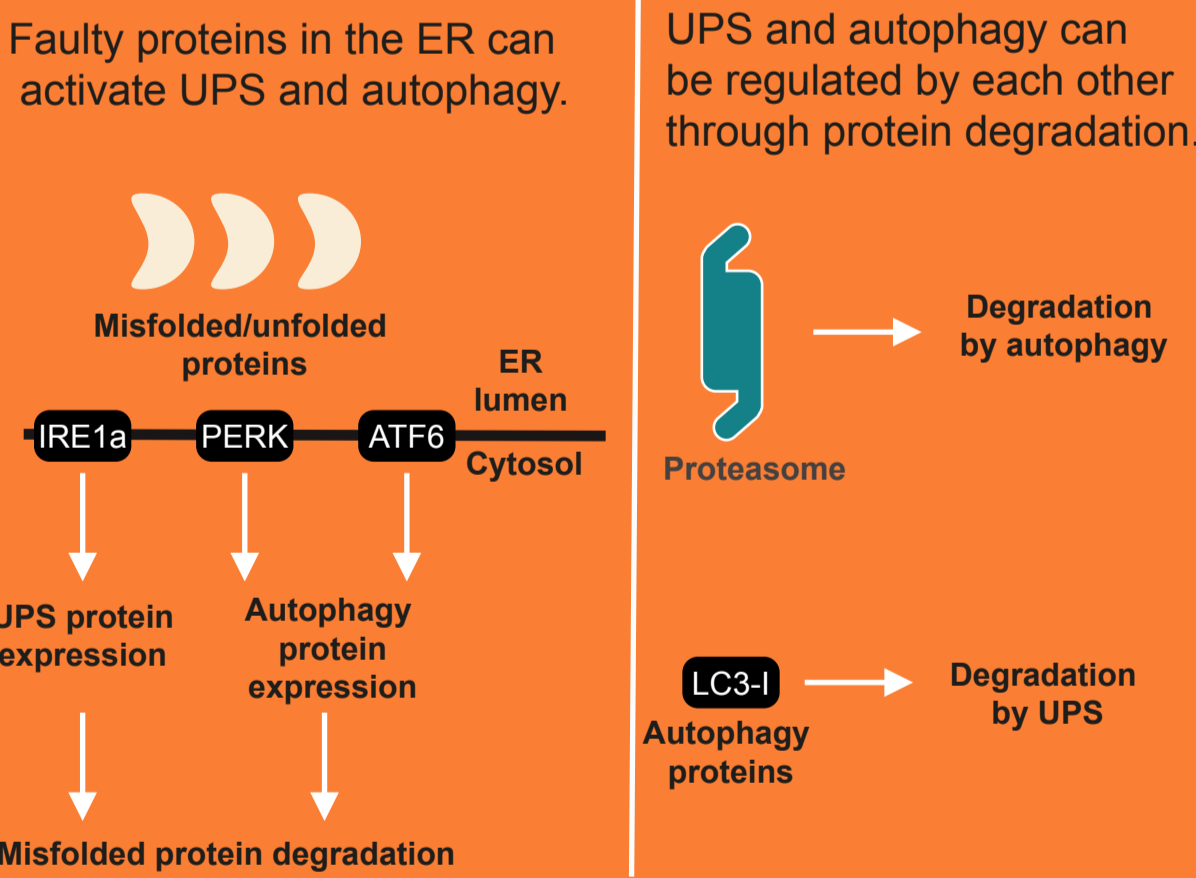
Ubiquitination of proteins, where proteins are modified with ubiquitin (Ub), is a process that relies on a series of enzymes: E1, E2, and E3 (such as MDM2). Multiple Ub molecules can be added to proteins, creating diverse ubiquitin 'tags' that determine the fate of proteins.



**Waste Processing Station:**  
Protein degradation



### Interconnecting Pathways



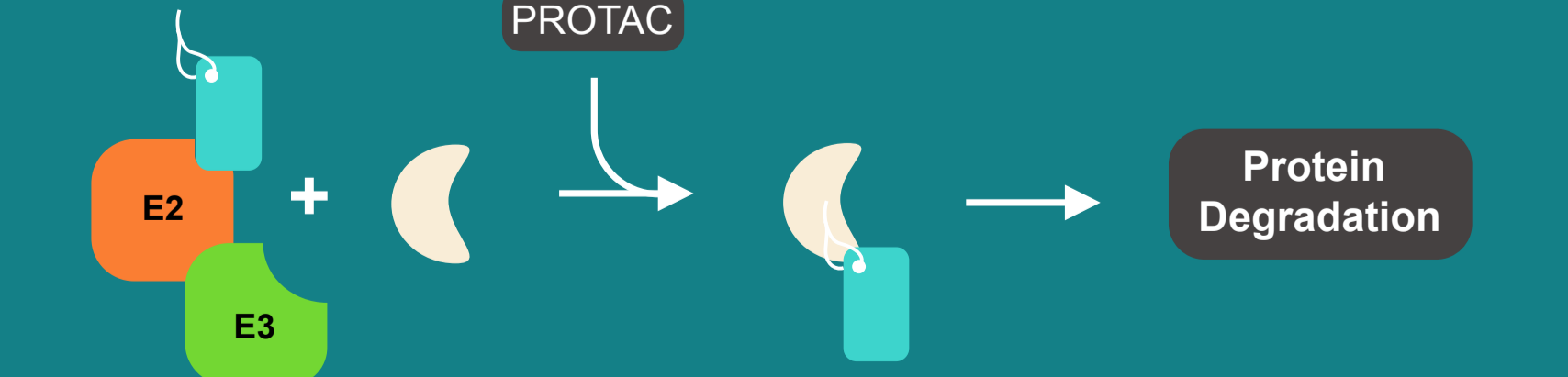
**Waste Collection:**  
Broken down protein products ready to be reused by cells

## Neurodegenerative disease and TPD

Misfolded proteins—such as Tau and TDP-43—can clump together, forming aggregates that are toxic to cells. If found in neurons, these aggregates can lead to neurodegenerative diseases such as Alzheimer's Disease and amyotrophic lateral sclerosis (ALS).



TPD involves hijacking protein degradation pathways to clear pathological proteins—which may provide a therapeutic strategy to treat neurodegenerative diseases. For example, TPD can use PROteolysis TArgeting Chimeras (PROTACs), which bind a target protein and E3 ligase. The target protein is ubiquitinated and degraded.



For more information on TPD, visit: [www.revivity.com](http://www.revivity.com)



### References

Nur Mehpare Kocaturk et al. Crosstalk Between Mammalian Autophagy and the Ubiquitin-Proteasome System. *Frontiers in Cell and Developmental Biology*. October 2018.  
 PerkinElmer. Protein Degradation Guide to the Proteasome and Autophagy Systems.  
 Yingxu Fang et al. Progress and Challenges in Targeted Protein Degradation for Neurodegenerative Disease Therapy. *Journal of Medicinal Chemistry*. August 2022.  
 Hwan-Ching Tai et al. Ubiquitin, the Proteasome and Protein Degradation in Neuronal Function and Dysfunction. *Nature Reviews Neuroscience*. December 2008.  
 Ze'ev A. Ronai. Monoubiquitination in Proteasomal Degradation. *PNAS*. August 2016.  
 Malene Hansen et al. Autophagy as a Promoter of Longevity: Insights from Model Organisms. *Nature Reviews Molecular Cell Biology*. July 2018.  
 M Catarina Silva et al. Targeted Degradation of Aberrant Tau in Frontotemporal Dementia Patient-Derived Neuronal Cell Models. *eLife*. March 2019.