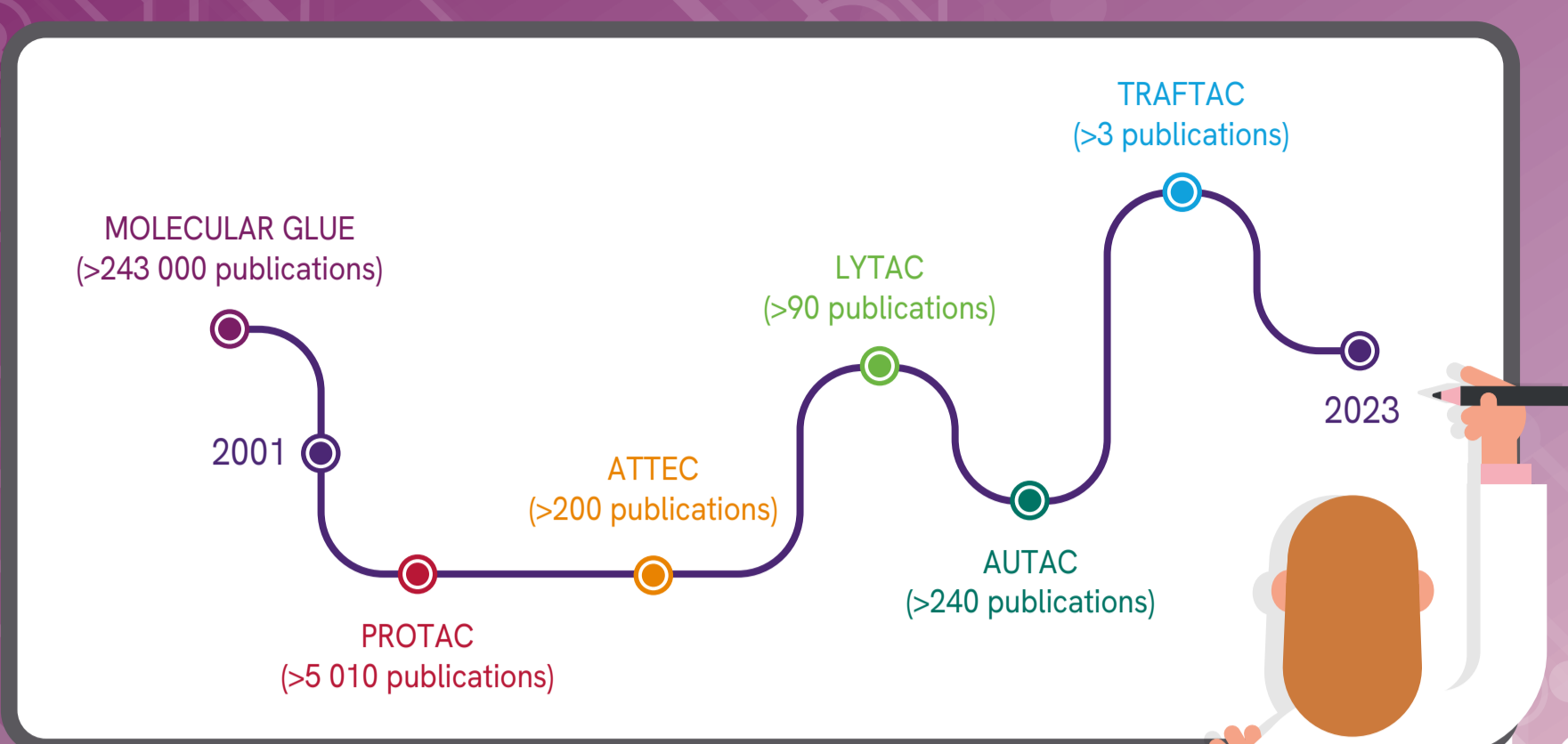


Targeted protein degradation (TDP), the next disruption in drug discovery?

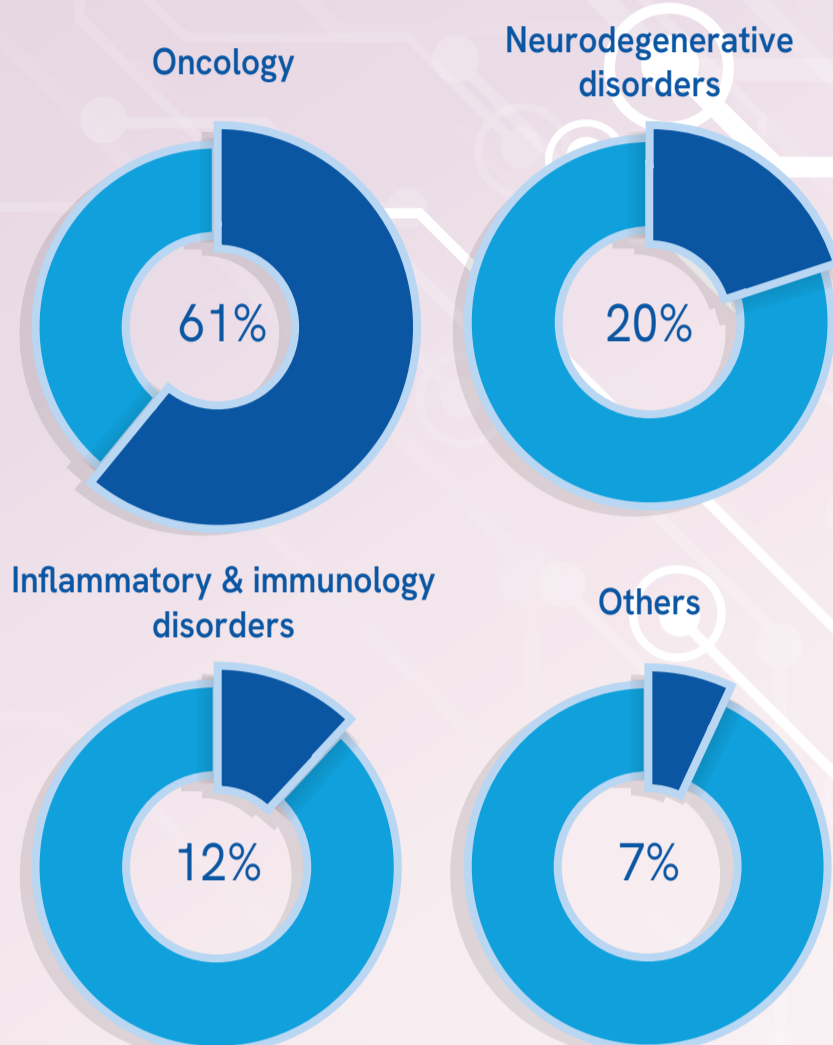


TDP technologies have been redefining the game for the past 20 years...



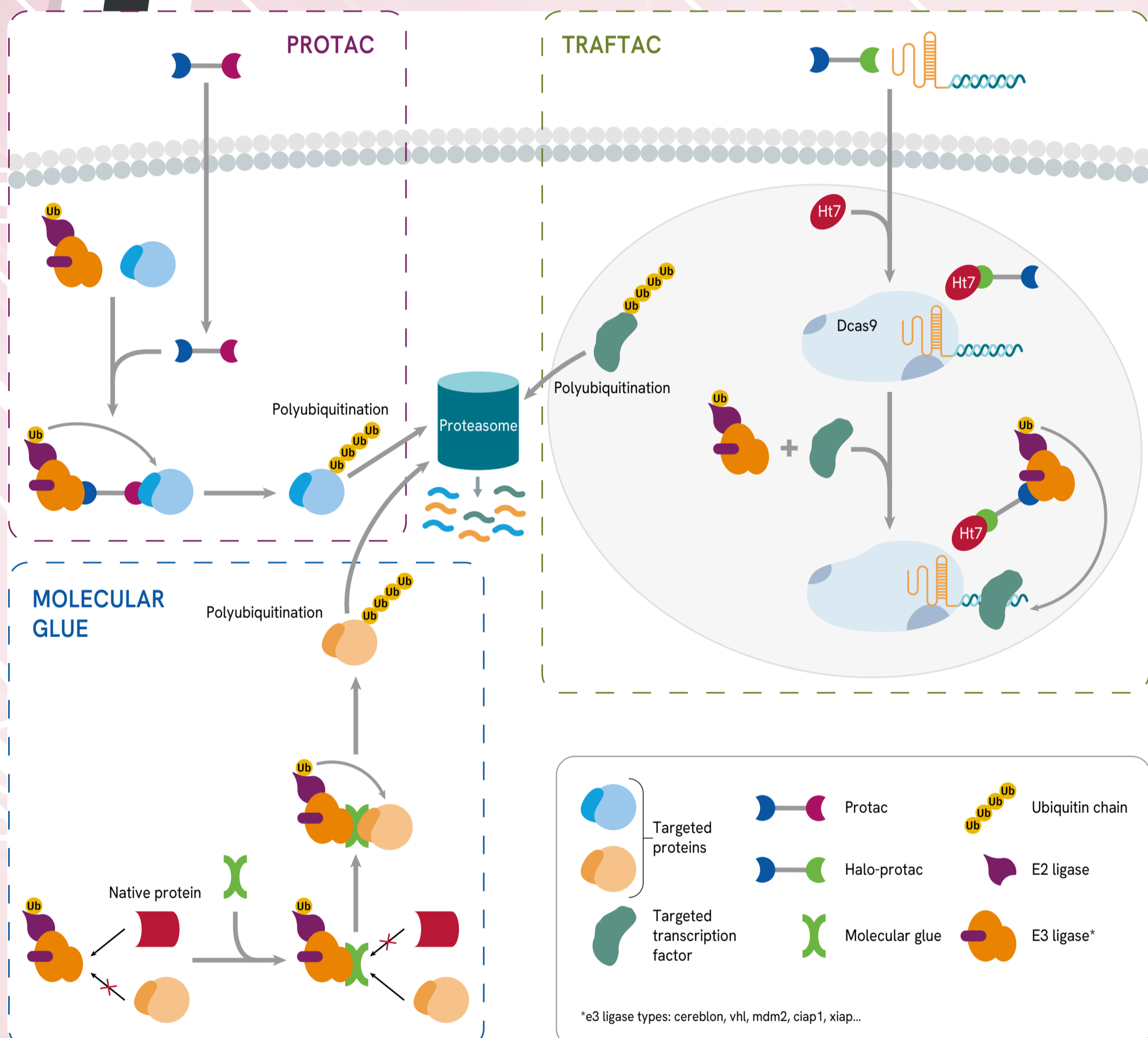
... they could cover a wide range of therapeutic areas...

... and solve new challenges.

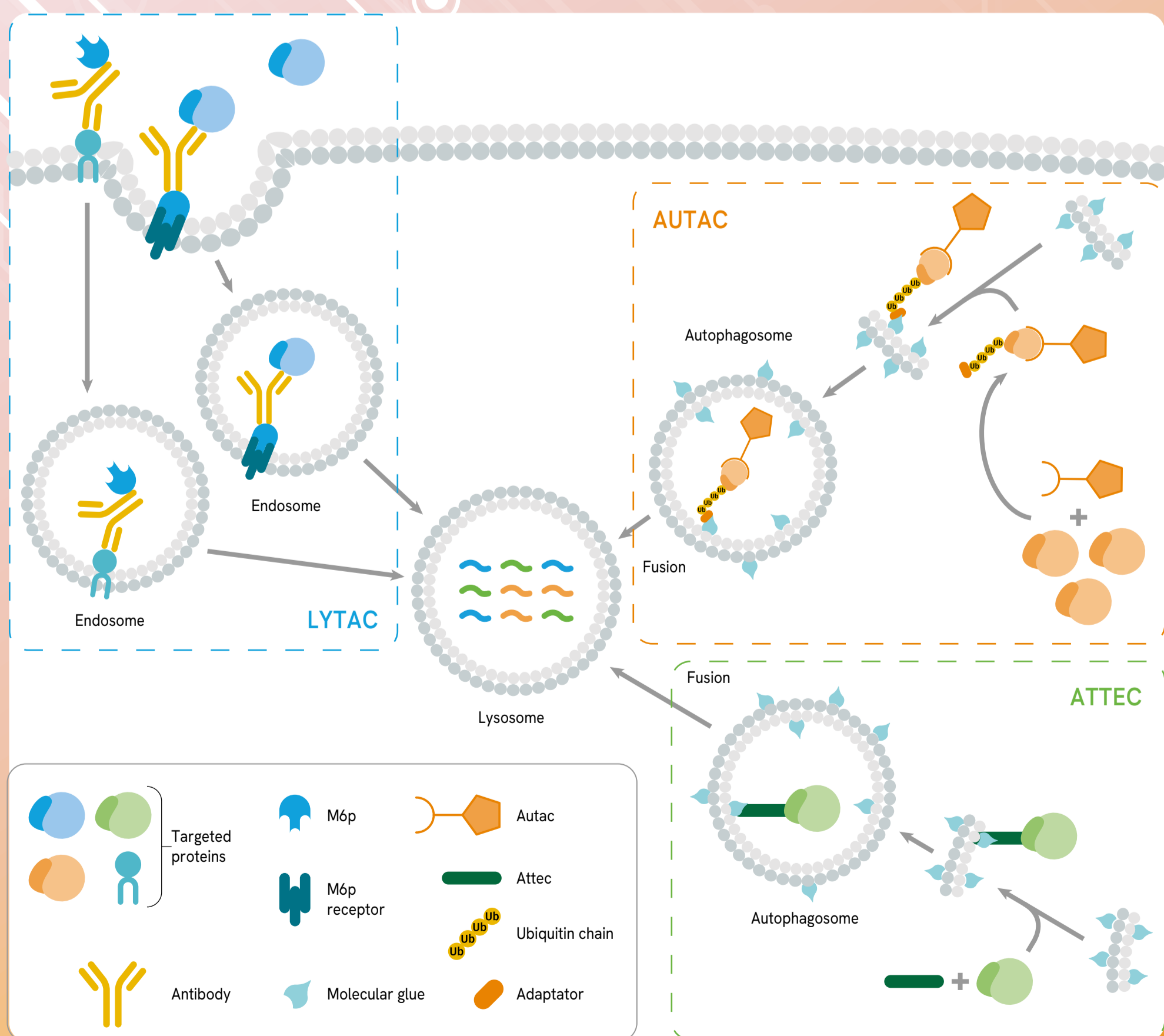


How do they work?

Ubiquitin proteasome system



Autophagy lysosomal system



References

- Yu ding, yihan fei, boxun lu. Emerging new concepts of degrader technologies. 2020.04.005
- Mikihiro naito, nobumichi ohoka, norihito shibata and yoshinori tsukumo. Targeted protein degradation by chimeric small molecules, protacs and snipers. Frontiers in chemistry. 2019.00849
- Momar toure, craig crews. Small-molecule protacs : new approaches to protein degradation. 201507978