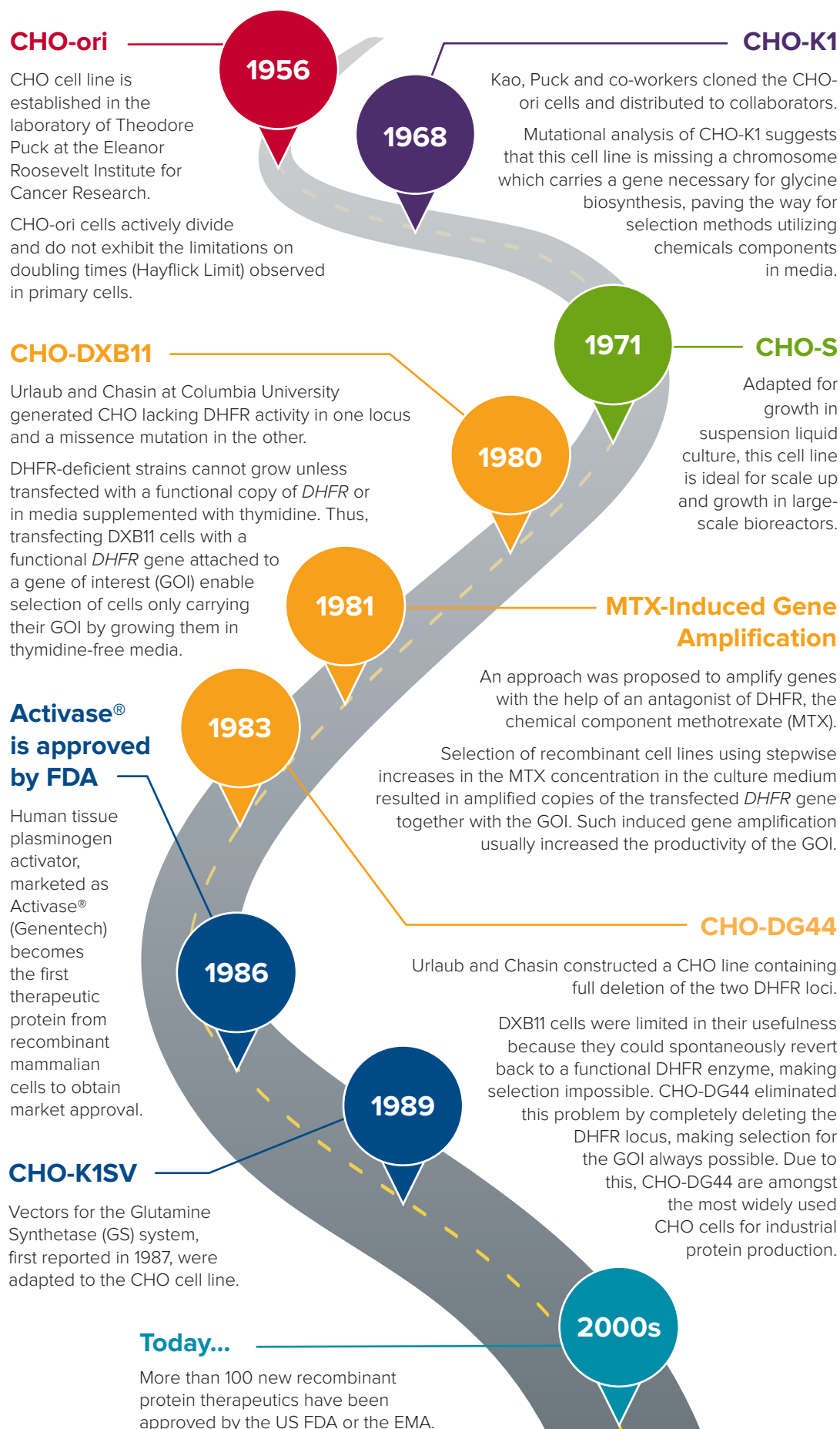




# The evolution of CHO cells' role in cell line development


Since the first approval of recombinant insulin and human growth hormone in the early 1980s, a multitude of recombinant protein therapeutics have been approved by regulatory agencies, notably the FDA in the US and EMA in Europe. Given this significant increase in successful introduction of biological therapeutic agents, there is a crucial need in the drug discovery space to support more efficient manufacturing processes that require highly productive cell lines. Consequently, **CHINESE HAMSTER OVARY (CHO)** cells have emerged as the gold standard for the manufacturing and regulatory approval of therapeutic proteins.

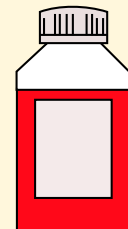
## MILESTONES

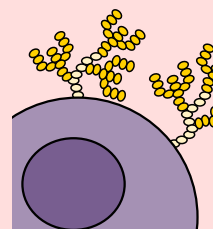


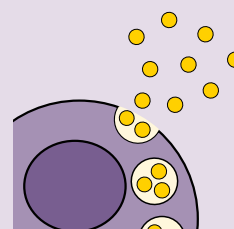
## WHY CHO?

Several key properties of CHO cells have driven their establishment as the preferred host cell line for regulatory approvals of recombinant therapeutic products:

- 

Adaptable to growing in suspension culture, which is ideal for large scale production in bioreactors.
- 

Adaptable to growth in serum-free and chemically defined (animal-free) media supplements, which ensures reproducibility between different batches of cell culture.
- 

Allow post-translational modifications (e.g. glycosylations) to recombinant proteins which are compatible and bioactive in humans.
- 

Several chemical selection and gene amplification systems have been developed for CHO cells, optimized for higher yield of recombinant protein per cell.

Cell line development can be time-consuming.

**ACCELERATE THE DISCOVERY WITH OUR PRODUCTS**



**CHO Growth A**  
Liquid and semi-solid culture media specifically for CHO cells



**ClonePix 2 System**  
Automated clone screening and picking system



**CloneSelect Imager**  
Imaging system to track and confirm monoclonality