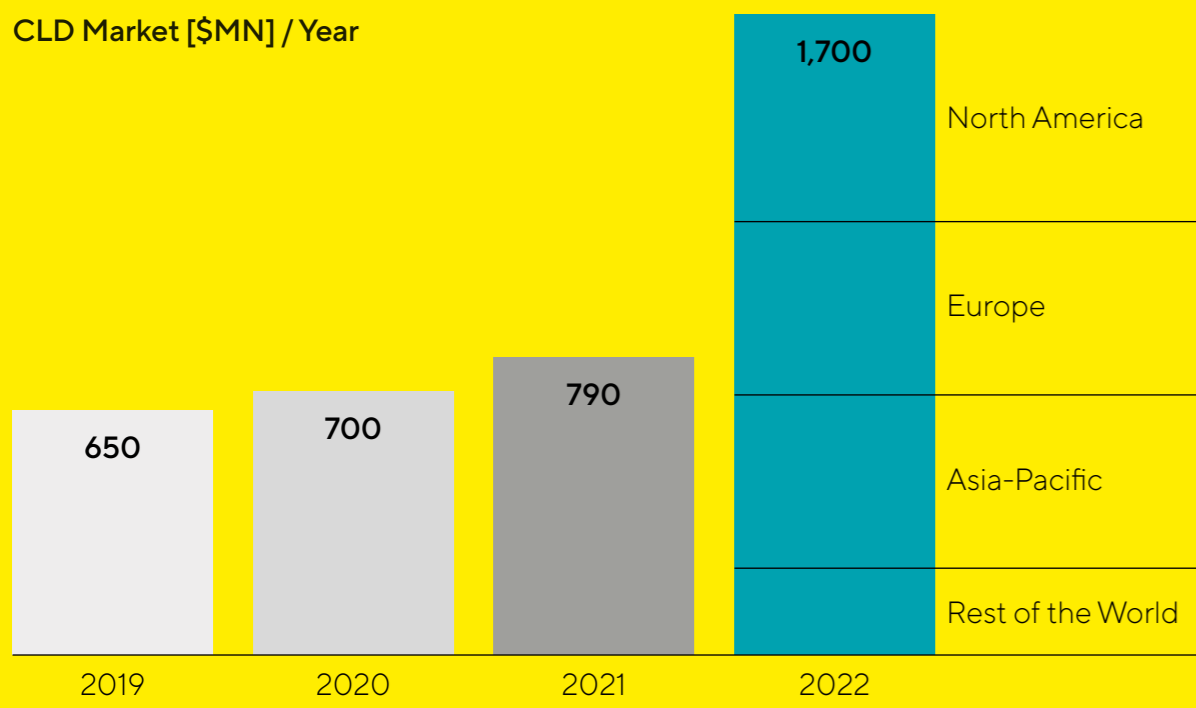


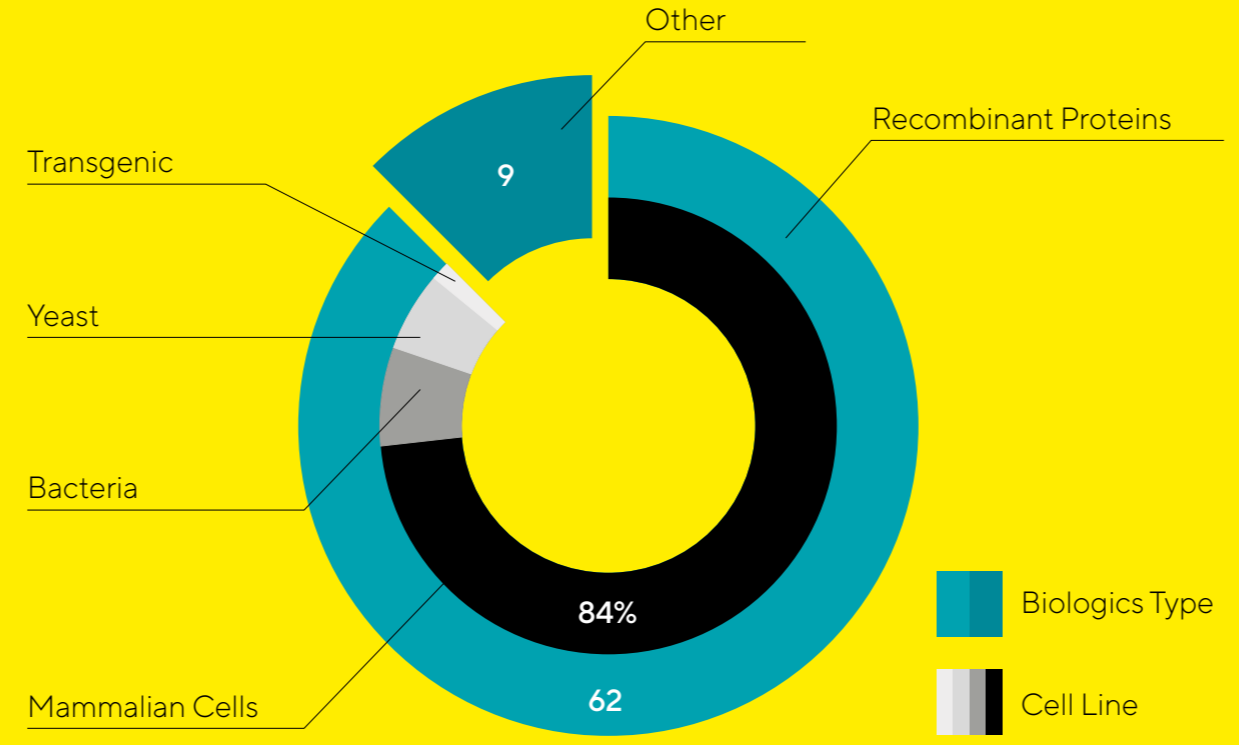
Cell Line Development Facts & Figures

The global cell line development service market is expected to reach **\$1.7 billion** in revenue by **2028**.

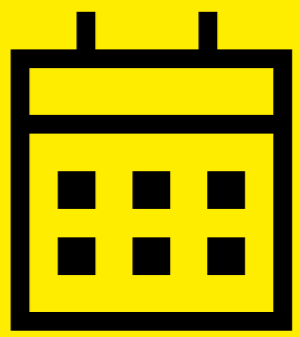
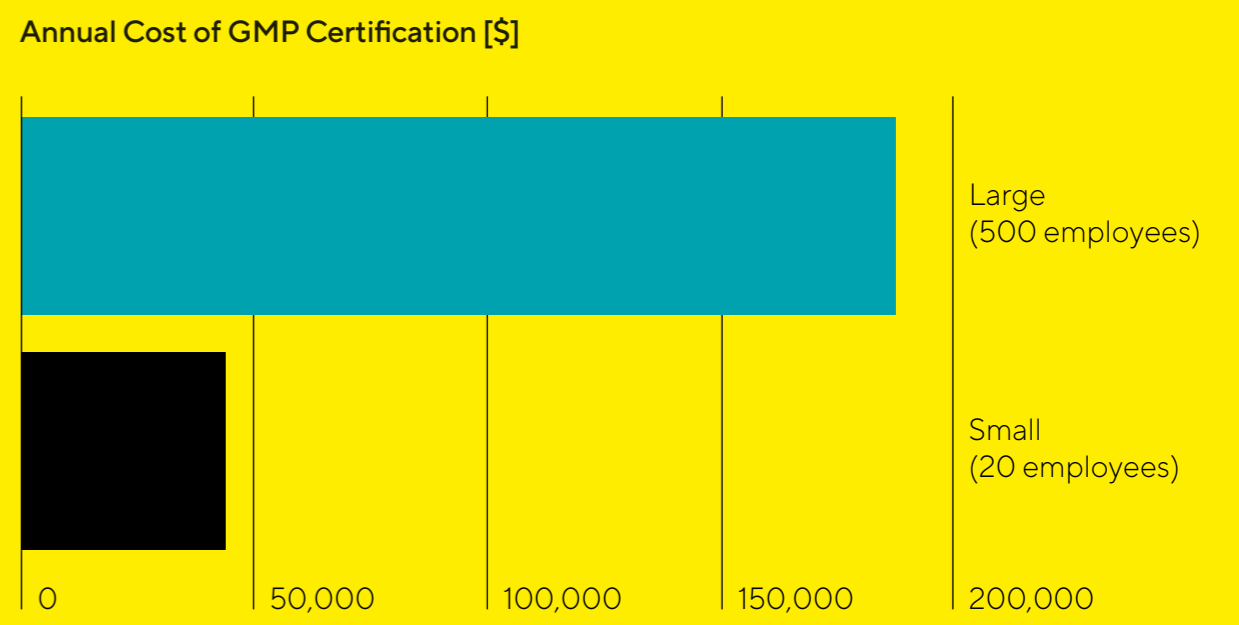


North America is expected to command over a **third** of the **market share**.

62 of the 71 newly approved biopharmaceutical active ingredients in the last 4 years were recombinant proteins and of those, **84%** were from **mammalian cells**.



GMP certification can be **pricey** – for small biotechs (<20 employees), the FDA estimates initial costs of around **\$26,000**, plus **\$46,000 annually**.



Cell line development is a **time-consuming** process – GMP grade cell lines are typically developed in **12-18 months**.

