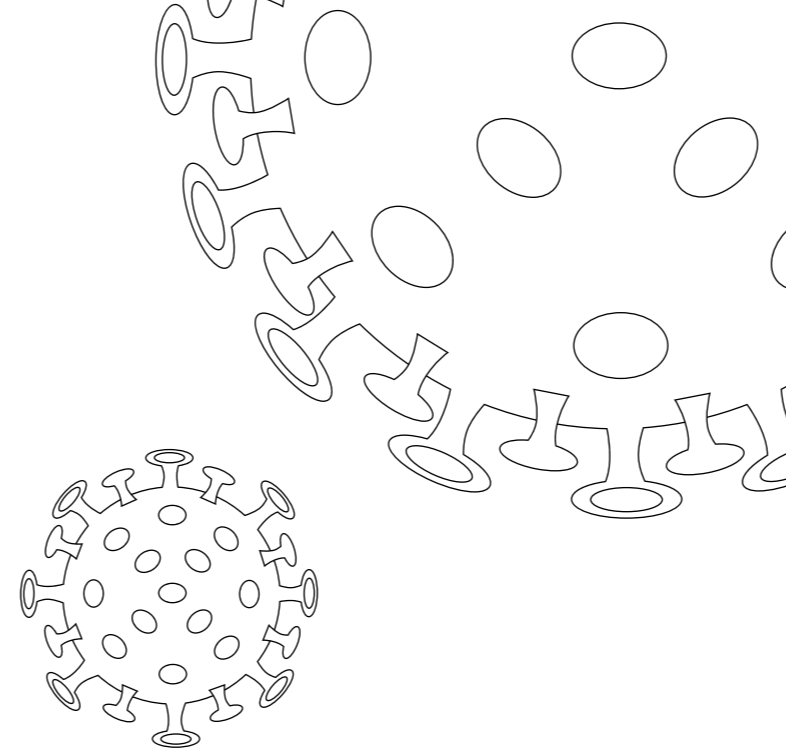


COVID-19 Survey on Learnings & Outlook for Future Pandemics

As the industry focuses its efforts to bring COVID-19 treatments swiftly to market, there is renewed interest in the development and manufacturing of vaccines. A survey has recently been conducted to determine what thoughts and opinions are regarding current and future vaccine development. Results of the survey are shown here.



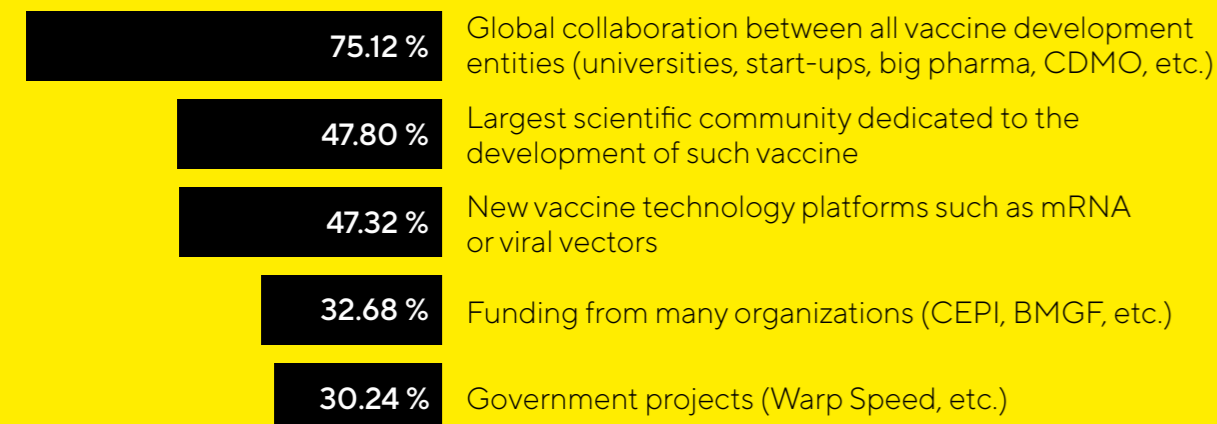
Do you think learnings and approaches from accelerated vaccine development of a vaccine against SARS-CoV-2 will be replicable to other vaccines needed for future outbreaks?

42.4%^{yes}

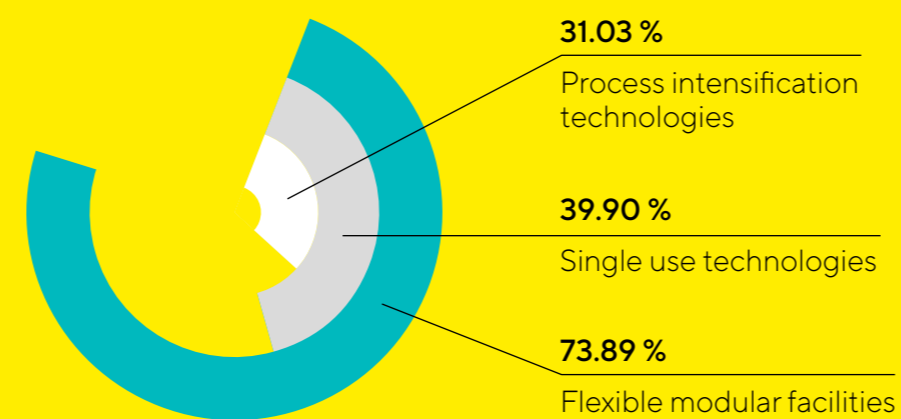
50.2%^{optimistic}

7.4%^{no}

What major factor(s) will allow to get a vaccine in less than 24 months? (select all that apply)



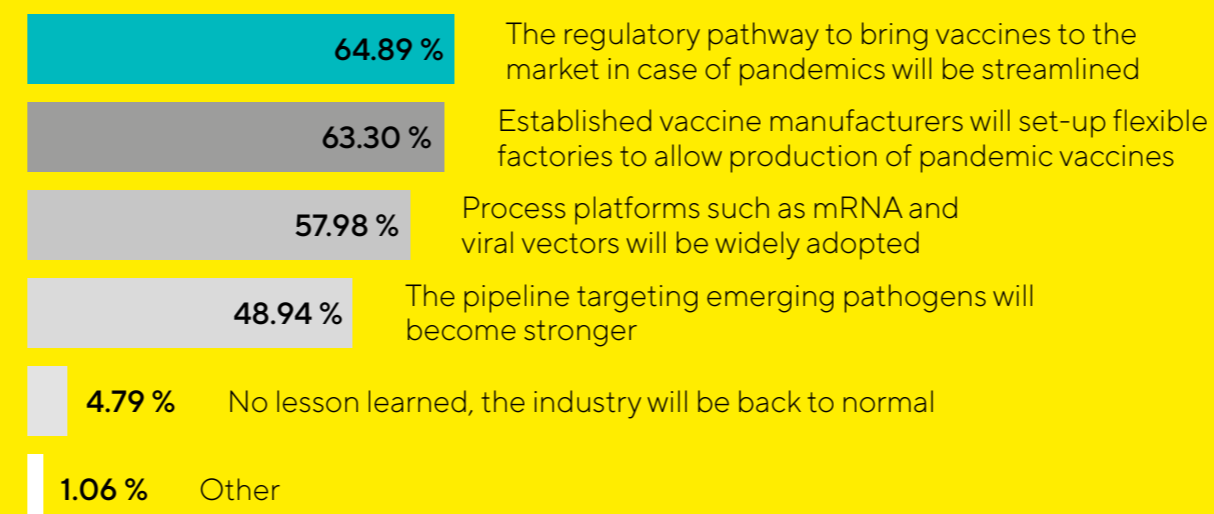
What breakthrough technology/service will allow building production capacity in the shortest timelines? (select all that apply)



To which extent will this pandemic change the vaccine developer/manufacturer behavior? (select all that apply)



How will the COVID-19 outbreak influence the future of the vaccine industry? (select all that apply)



What is the main challenge in bringing the vaccine to market under severe time pressure?

