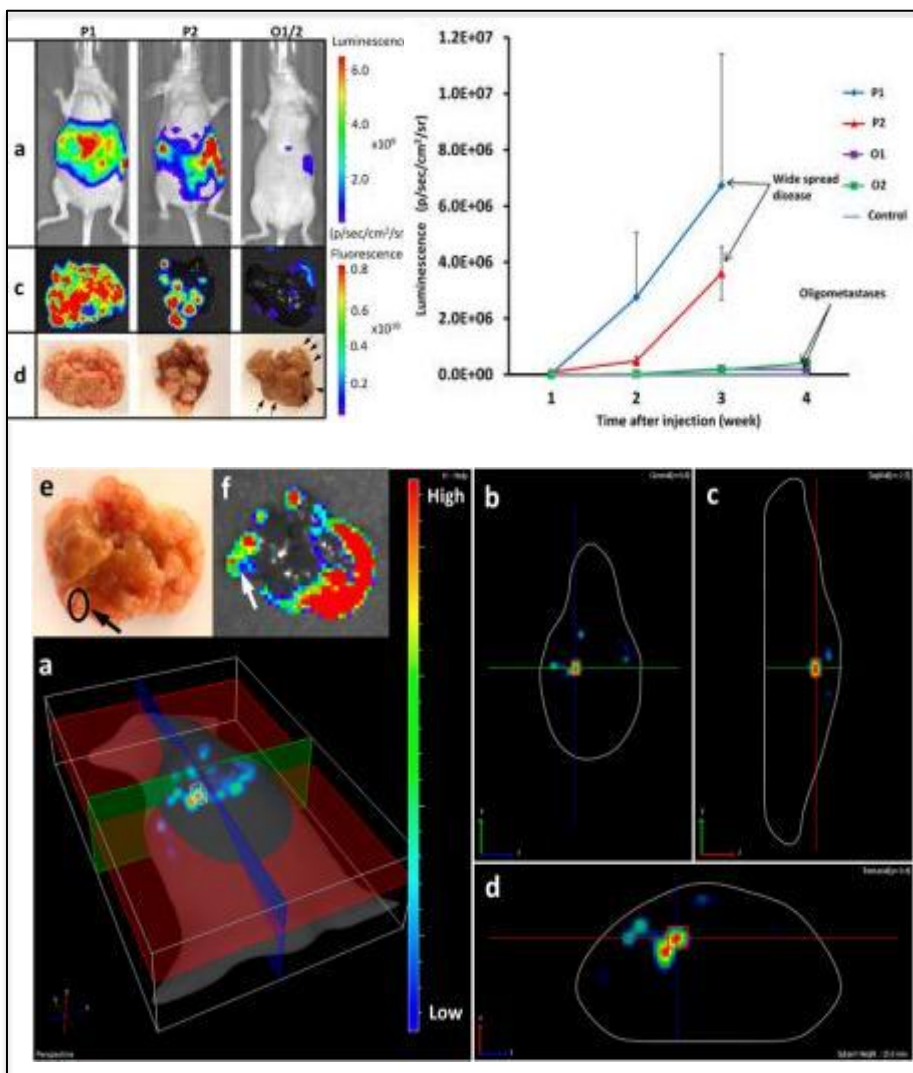
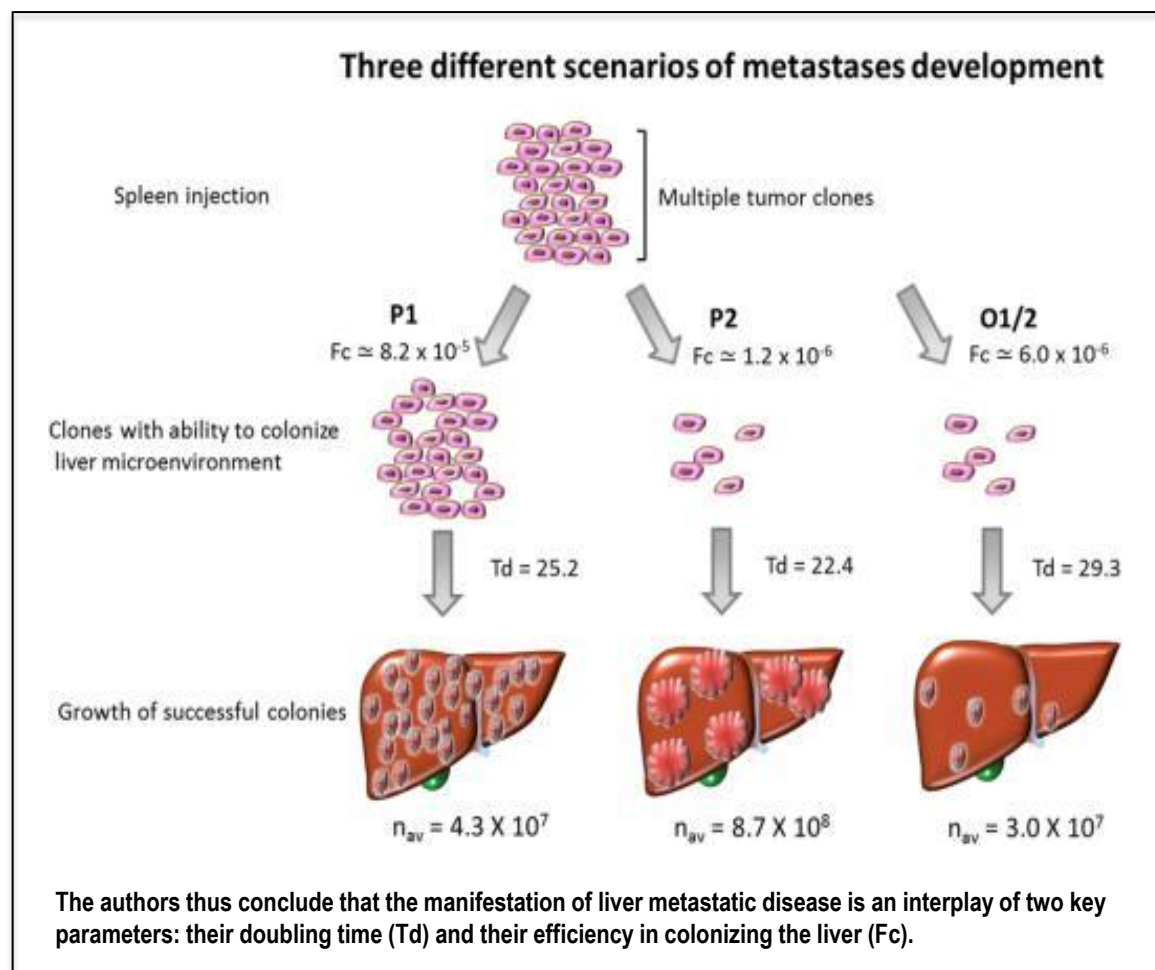


Colorectal cancer patients often develop liver metastases and thus, frequently have poor prognosis. There additionally appears to be a vast heterogeneity in their liver metastatic disease, a characteristic that hasn't been adequately explored in animal models of colorectal cancer. While bioluminescence imaging (BLI) has been widely used to non-invasively monitor colorectal cancer and liver metastatic development *in vivo*, a study specifically emphasizing their growth rates and colonization efficiencies within the liver microenvironment hasn't been attempted until now.



Imaging was used to screen 16 individual clones and identify 2 polymetastatic clones (P1 and P2) that form widely disseminated metastatic tumors and 2 oligometastatic clones (O1/2) characterized by a limited metastatic phenotype. Further quantification of their growth and colonization ability by imaging revealed:

- P1 has a high colonization potential but a low growth rate → large number of small metastatic tumors
- P2 has lower colonization capacity but faster doubling time → small number of large secondary tumors
- O1/2 show lower colonization efficiencies and lower growth → limited metastatic disease (oligometastatic)



The authors used a combination of dual optical reporters (luciferase and tdTomato), *in vivo* 2D BLI and *ex vivo* analysis to longitudinally detect, quantify and resolve metastatic colonies. 3D bioluminescence tomography (DLIT) was used to further validate the spatiotemporal dynamics of metastatic development.