

➤ SubQperior and orthotopic breast cancer model

JIMT1 origin is a breast tumor. Therefore subQperior is a special case: it is orthotopic. The method for tumor cell implantation into the mammary fat pad is the same procedure as orthotopic implantation of breast cancer cells. JIMT1 cells do not express luciferase.

➤ Jimt1 cells

Human JIMT1 cells were isolated from a patient with a invasive ductal carcinoma.

As routine quality controls, the cells are regularly checked for Mycoplasma contamination and authenticity (via STR DNA Typing).

➤ Tumor growth *in vivo*

JIMT1 cells harvested from tissue culture flasks are implanted into the mammary fat pad of the mice. Resulting tumors are monitored by calipering twice weekly.

Animal weights are measured three times weekly.

Animal behaviour is monitored daily.

All mice are maintained in separated isolated housing at constant temperature and humidity.

Accessory services: tumor wet weight and volume measurement at necropsy, blood sampling, paraffin embedding of tumor tissue, histological & pathological analysis, provision of tumor tissue for target validation.

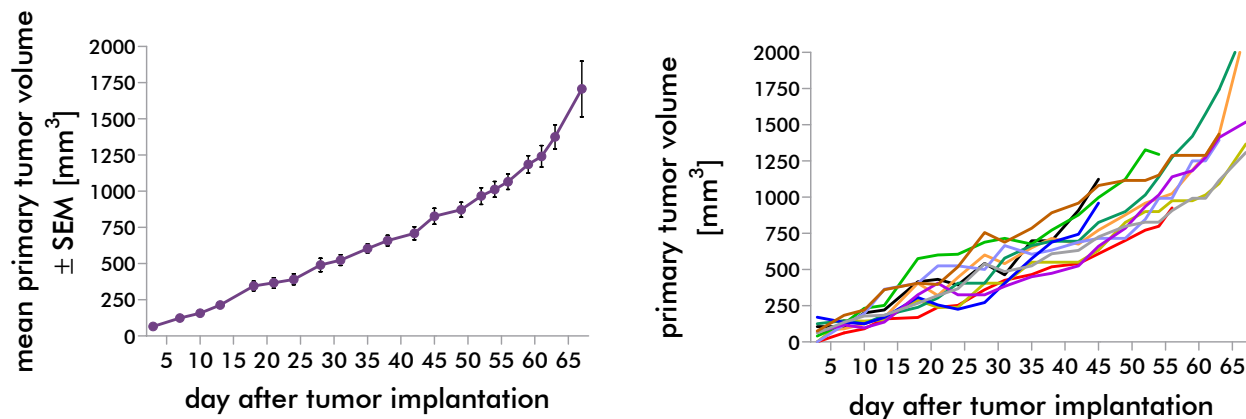


Figure 1: Tumor growth of JIMT1 cells in a subQperior xenograft in vivo. Left graph: mean values. Right graph: individual tumors.

➤ Study example

If you are interested in receiving information on potential positive controls please reach out to our Business Development team at requests@reactionbiology.com.