

# SubQperior

the next generation of tumor models

A superior implantation method for cell-line derived tumor models

- Homogeneous tumor growth
- Reproducible study outcome
- Outstanding statistical value

Standard implantation via subcutaneous injection causes frequent ulceration resulting in early abrogation of studies, leaving researchers with too short treatment windows and high heterogeneity with poor statistical value of study results.

To overcome these challenges, we have developed a superior implantation method for standard cell-line derived tumor models.

subQperior: tumor cell implantation into the mammary fat pad.

A decorative graphic consisting of seven light purple circles of varying sizes arranged in a grid-like pattern on the right side of the slide.

Let's discover together.









