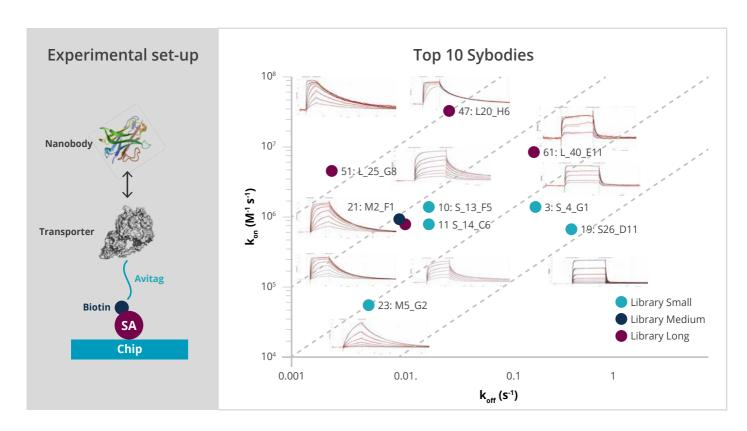
SYBODIES CAN SHOW GREAT DIVERSITY IN KINETIC PROPERTIES





"The combination of unique technology, robustness, flexibility in designing a broad range of assays and high sensitivity, enabling even fragment sized molecules to be screened and characterized, made the Creoptix WAVEdelta the instrument investment of our choice. Thorough testing in our laboratory has convinced us that Creoptix' WAVE technology aids us best in discovering better drug molecules"

Nicolas Bocquet, Senior Scientist at



