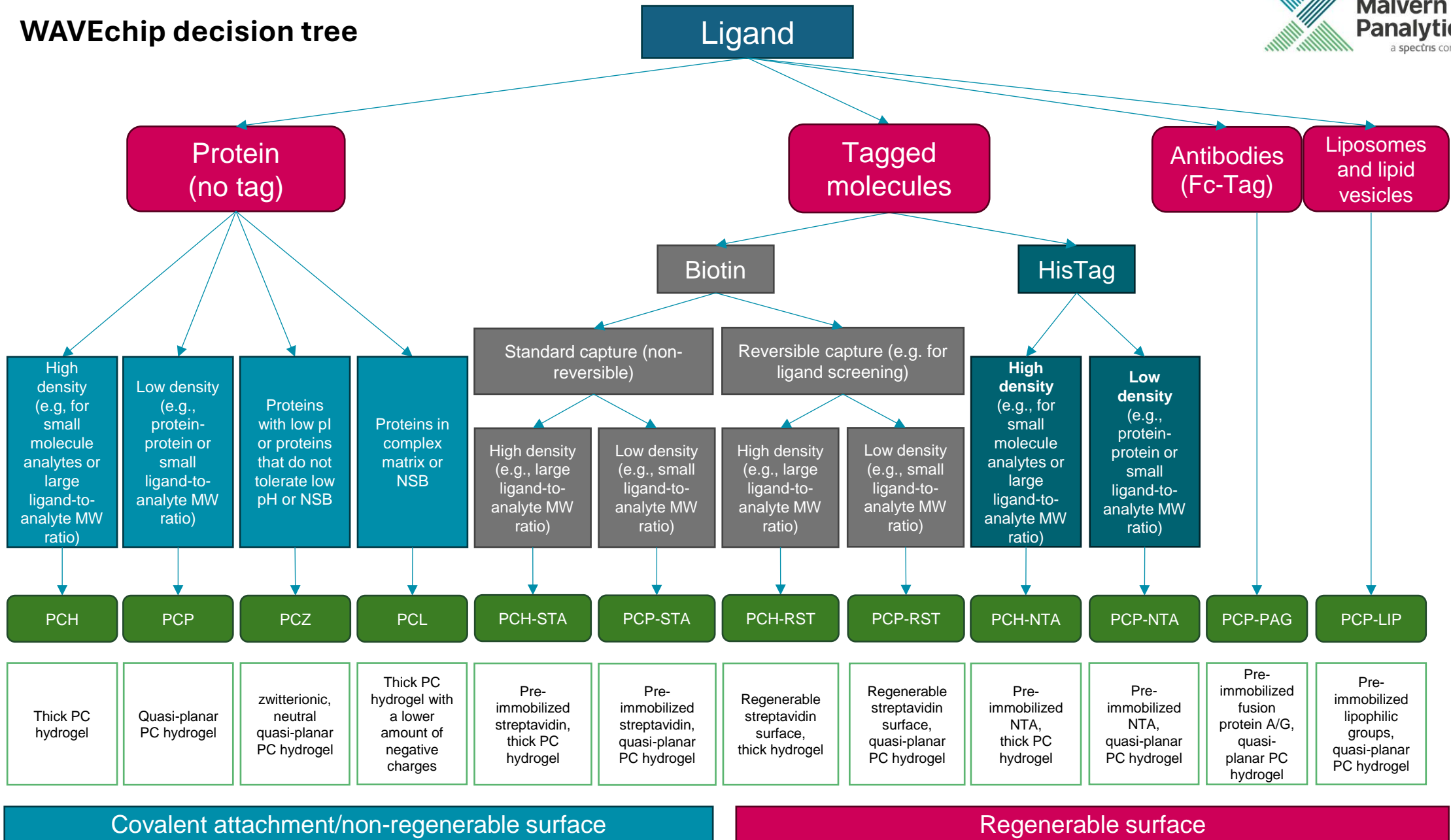


WAVEchip decision tree



Covalent attachment/non-regenerable surface

Regenerable surface