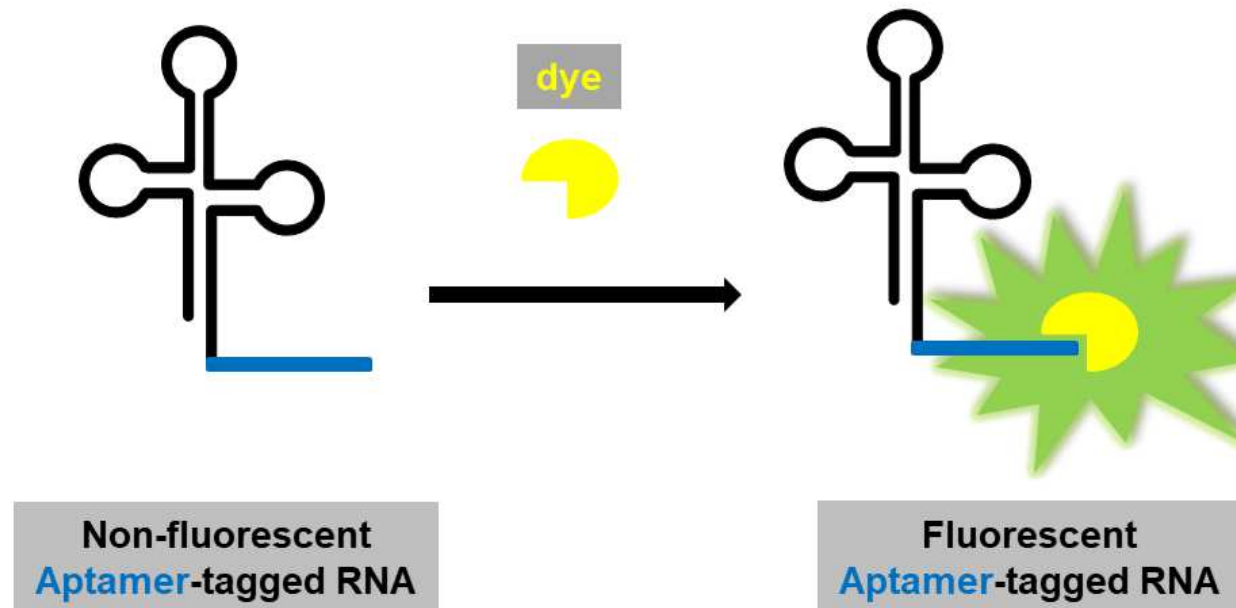


## Principle of light-up aptamer-based RNA imaging & analysis



**Figure 1** A specific aptamer sequence is attached to the 5'-, 3'- or internal position of a RNA sequence of interest e.g. by *in vitro* transcription. Once an aptamer specific dye is added, a fluorescence signal can be detected due to dye activation upon non-covalent binding to its corresponding aptamer sequence