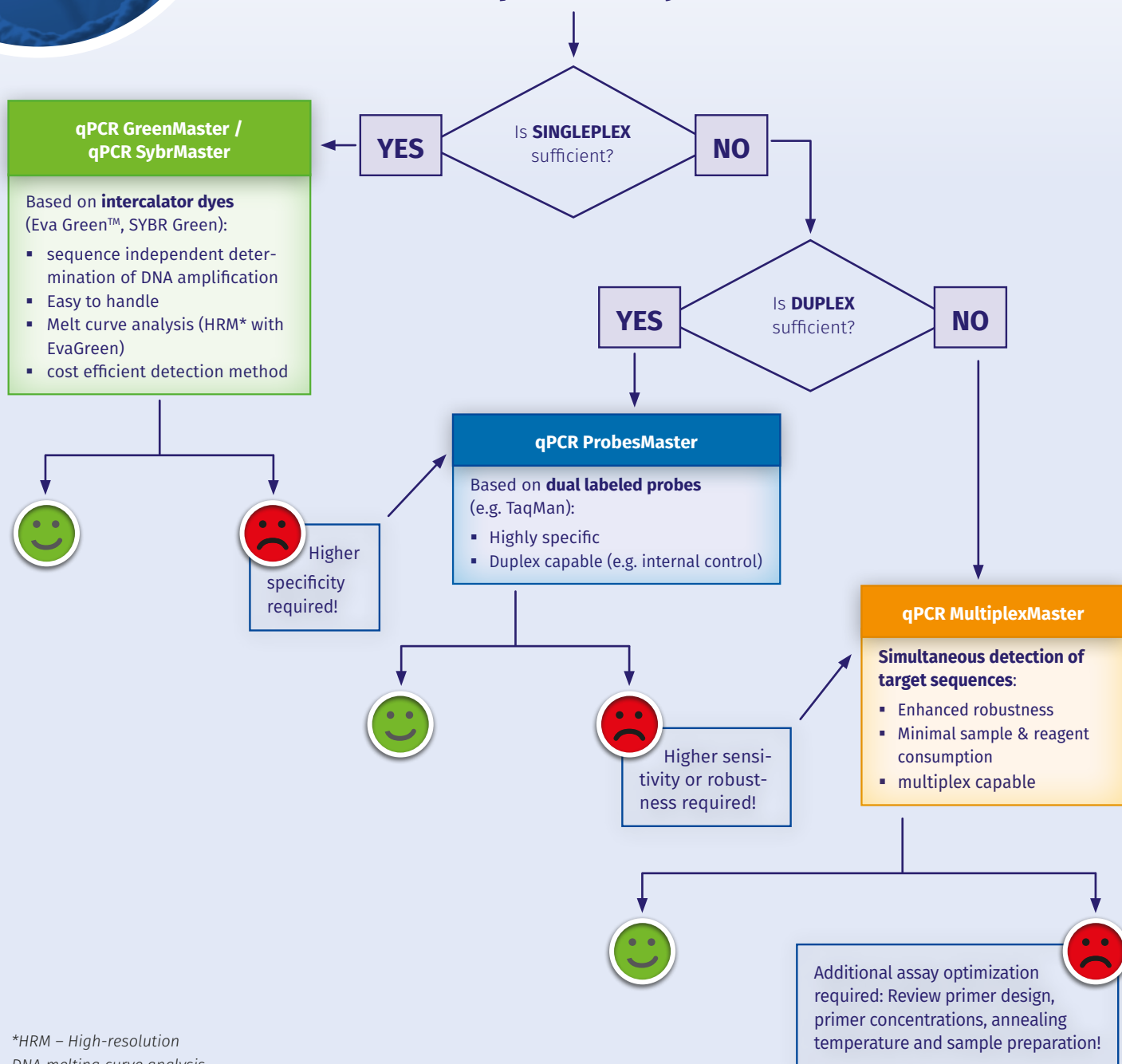


qPCR-mixes: Which one to use?



Quantitative Real-Time PCR (qPCR) is commonly applied to assess nucleic acid quantities. Its design & set up, optimization and potential trouble-shooting however, can be time-consuming, expensive and simply unnecessary! Learn - in this short tutorial - how to set-up the most promising assay.

Sample to Analyze



*HRM – High-resolution DNA melting curve analysis

