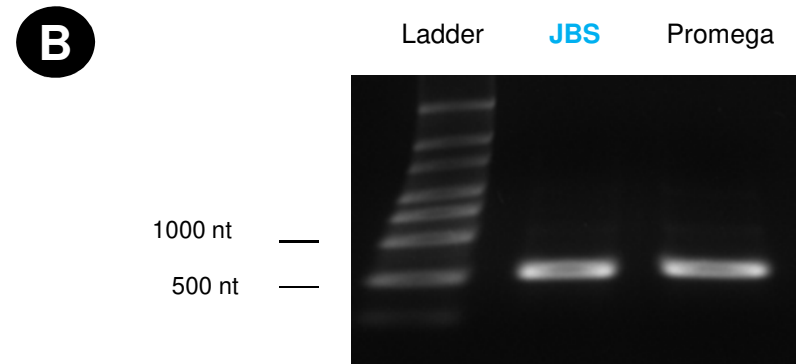
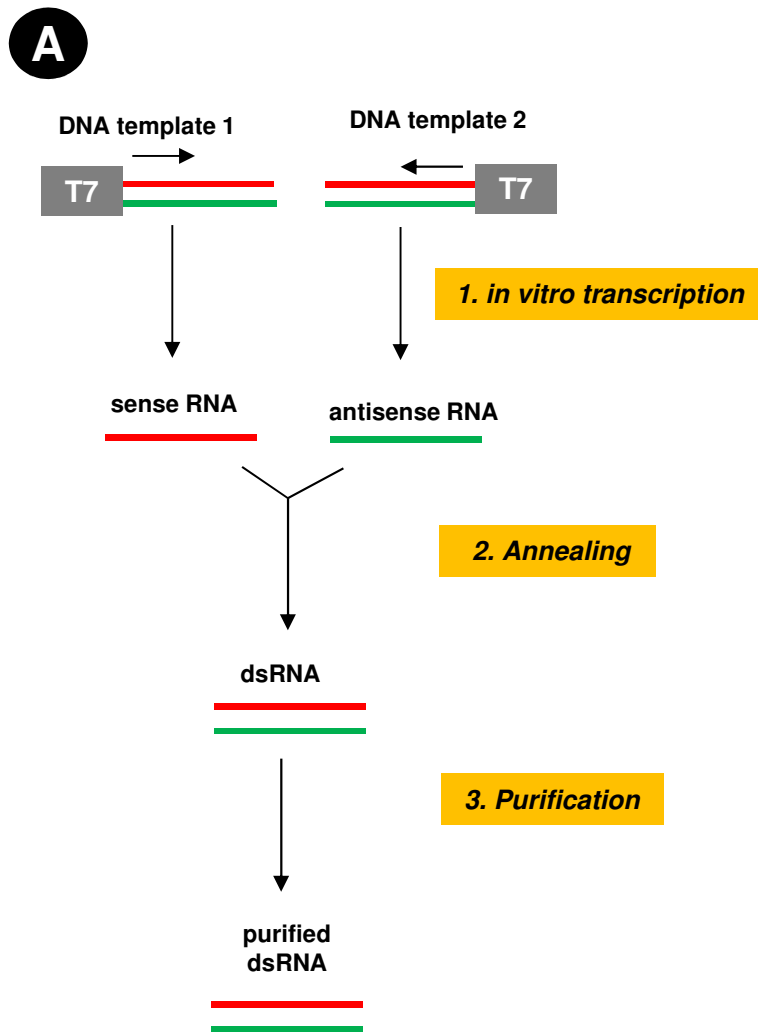


Similar results with Jena Bioscience's HighYield T7 RNAi Kit and the T7 RiboMAX™ Express RNAi System



C

Product	Cat. No.	Amount	Price €	Average yield / 20 µl reaction
HighYield T7 RNAi Kit	RNT-134	50 reactions à 20 µl	265	> 50 µg (> 2 mg / ml reaction)
T7 RiboMAX™ Express RNAi System	P1700	50 reactions à 20 µl	547	> 50 µg (> 2 mg / ml reaction)

Figure 1 Similar results with Jena Bioscience's HighYield T7 RNAi Kit and the T7 RiboMAX™ Express RNAi System.

A) dsRNA is synthesized in a 3 step procedure.

B) Similar transcription efficiency of Jena Bioscience's HighYield T7 RNAi Kit and T7 RiboMAX™ Express RNAi System. *In vitro* transcription has been performed according to the manufacturer protocol using 1 µg DNA template 1. Transcription efficiency was analyzed by agarose gelelectrophoresis (equal volumes of *in vitro* transcription reactions). JBS: HighYield T7 RNAi Kit, Promega: T7 RiboMAX™ Express RNAi System

C) Competitive pricing of Jena Bioscience's HighYield T7 RNAi Kit.