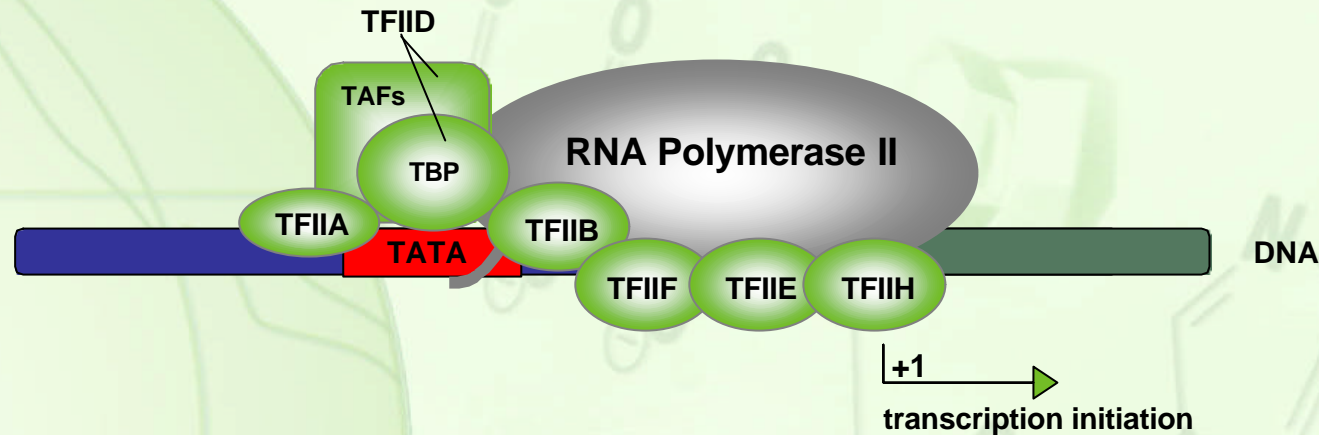


Transcription pre-initiation complex – basal factors of transcription by RNA polymerase II



The various transcription factors II and/or their subunits are available from Jena Bioscience!

- The eukaryotic pol II promoter consists of core elements which include the TATA box and other DNA sequences that define transcription start sites, and regulatory elements which either enhance or repress transcription in a gene-specific manner
- The core promoter is the site for assembly of the transcription pre-initiation complex which includes RNA pol II and the general transcription factors TBP, TFIIA, TFIIB, TFIIE, TFIIF, and TFIIH