

Tagmentase

Cat. No. C01070010

Lot: 2/a

Format: 10 μ l / 20 μ l

Concentration: 2 mg/ml

Product description

Diagenode Tagmentase is a hyperactive transposase with the ability to cut DNA and insert sequences of interest into any target DNA in one step. This enzyme is not loaded with DNA oligos.

Storage conditions

Store at -20°C . Guaranteed stable for 6 months from date of receipt when stored properly.

Storage buffer

Supplied in solution containing 50% v/v glycerol.

Precautions:

This product is for research use only. Not for use in diagnostic or therapeutic procedures.

Quality control

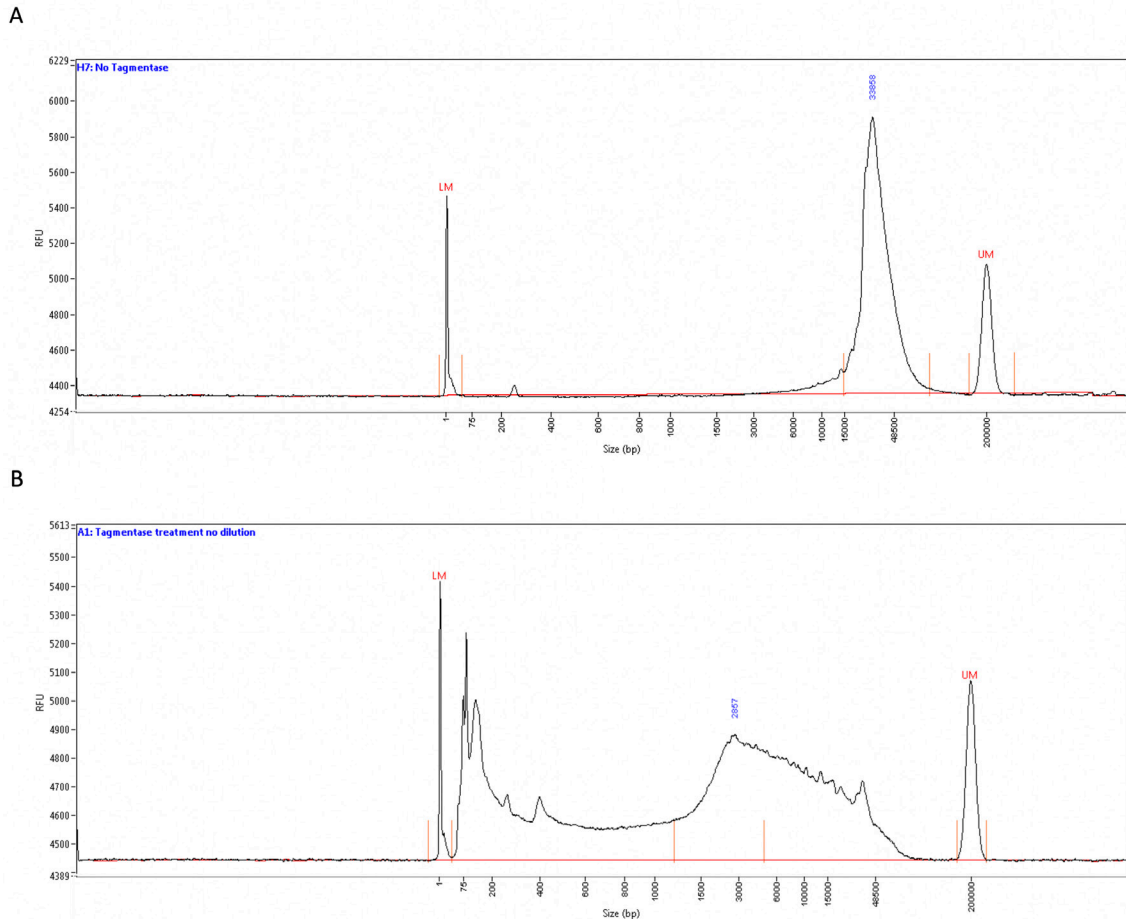


Figure 1: Efficient fragmentation of the lambda DNA after incubation with the Tagmentase

For fragmentation, 100 ng of DNA from bacteriophage lambda were incubated with diluted Diagenode Tagmentase (Cat. No. C01070010) for 7 min at 55°C. The reaction was stopped by addition of SDS (0.2% final concentration). After clean-up using AMPure XP beads (Beckman Coulter) on Diagenode IP-Star Compact Automated System, the size of the DNA was assessed on Fragment Analyzer (Agilent), using the HS Large Fragment 50kb Kit (Agilent). Profiles show the size of lambda DNA before (A) and after treatment with Tagmentase (B).



Figure 2: Fragmentation efficiency depending on the amount of Tagmentase

For fragmentation, 100 ng of DNA from bacteriophage lambda were incubated with Diagenode Tagmentase (Cat. No. C01070010) for 7 min at 55°C. The Tagmentase was previously diluted with the Tagmentase Dilution Buffer (Cat. No. C01070011) at 1/4 and 1/16 dilutions. The reaction was stopped by addition of SDS (0.2% final concentration). After clean-up using AMPure XP beads (Beckman Coulter) on Diagenode IP-Star Compact Automated System, the size of the DNA was assessed on Fragment Analyzer (Agilent), using the HS Large Fragment 50kb Kit (Agilent). The migration of the samples shows variations of the size distribution according to the amount of Tagmentase used for the reaction.