

Kpn I



Source: *Klebsiella pneumoniae* OK8.

Cat.-No.	Size	Conc.
EN-119S	3,500 units	10 u/μl
EN-119L	17,500 units	10 u/μl

Buffer supplied: 10x *Kpn* I and 10x BSA.

Substrate for unit definition:

λ DNA, *EcoR* I digest (2 sites).

Reaction conditions:

10 mM Tris-HCl (pH 7.0), 10 mM MgCl₂, 1 mM dithiothreitol, 0.01% Triton X-100, 100 μg/ml BSA. Incubate at 37°C.

Storage buffer:

50 mM KCl, 10 mM Tris-HCl (pH 7.4), 0.1 mM EDTA, 1 mM dithiothreitol, 200 μg/ml BSA and 50% glycerol.
Store at -20°C.

Ligation and recutting:

After 10-fold overdigestion with *Kpn* I, >95% of the DNA fragments can be ligated and recut with this enzyme.

Star activity:

Conditions of low ionic strength, high enzyme concentration, glycerol concentration >5% or pH >8.0 may result in star activity.

Heat inactivation: No.