

Alu I



Source: *Arthrobacter luteus*.

Cat.-No.	Size	Conc.
EN-101S	600 units	10 u/μl
EN-101L	3,000 units	10 u/μl

NB:
BSA is already included in the Buffer solution!

Buffer supplied: 10x B1 (incl. 10x BSA).

Substrate for unit definition: λ DNA (143 sites).

Reaction conditions:

10 mM Tris-HCl (pH 7.9), 10 mM MgCl₂, 1 mM dithiothreitol, 100 μg/ml BSA.

Incubate at 37°C.

Storage buffer:

100 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 *Alu I*, >95% of the DNA fragments can be ligated and recut with this enzyme.

Heat inactivation: 65°C for 20 minutes.