

SseB I

(Stu I)



Source: *Streptomyces* species.

Cat.-No.	Size	Conc.
EN-138S	1,500 units	10 u/μl
EN-138L	7,500 units	10 u/μl

NB:
BSA is already included in the Buffer solution!

Buffer supplied: 10x B3 (incl.10x BSA).

Substrate for unit definition:
λ DNA, *Hind* III digest (6 sites).

Reaction conditions:
100 mM NaCl, 50 mM Tris-HCl (pH 7.9), 10 mM MgCl₂,
1 mM dithiothreitol, 100 μg/ml BSA.
Incubate at **37°C**.

Storage buffer:
50 mM KCl, 10 mM Tris-HCl (pH 7.9), 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 SseB I, >95% of the
DNA fragments can be ligated and recut with this
enzyme.

Heat inactivation: 65°C for 20 minutes.