

Sph I



Source: *Streptomyces phaeochromogenes*.

Cat.-No.	Size	Conc.
EN-137S	250 units	10 u/μl
EN-137L	1,250 units	10 u/μl

NB:
BSA is already included in the Buffer solution!

Buffer supplied: 10x B2 (incl. 10x BSA).

Substrate for unit definition: λ DNA (1 site).

Reaction conditions:

50 mM NaCl, 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, 400 μg/ml BSA, and 50% glycerol.
Store at -20°C.

Ligation and recutting:

After 10-fold overdigestion with Sph I, >98% of the DNA fragments can be ligated and recut with this enzyme.

Heat inactivation: 65°C for 20 minutes.