

BstE II 5'...GGTNACC...3'
3'...CCANTGG...5'

Source: *Bacillus stearothermophilus*.

Cat.-No.	Size	Conc.
EN-144S	1,750 units	10 u/μl
EN-144L	8,750 units	10 u/μl

Buffer supplied: 10x *BstE II* and 10x BSA.

Substrate for unit definition: λ DNA (13 sites).

Reaction conditions:

100 mM KCl, 10 mM Tris-HCl (pH 7.4), 5 mM MgCl₂, 1 mM dithiothreitol, 0.1% Triton X-100, 100 μg/ml BSA.

Incubate at **60°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 40-fold overdigestion with *BstE II*, >95% of the DNA fragments can be ligated and recut with this enzyme.

Note:

BstE II exhibits 10-15% activity at 37°C.

Heat inactivation: No.