

***Bsi*S I**
(*Hpa* II)



Source: *Bacillus stearothermophilus*.

Cat.-No.	Size	Conc.
EN-111S	2,200 units	10 u/μl
EN-111L	11,000 units	10 u/μl

Buffer supplied: 10x *Bsi*S I and 10x BSA.

Substrate for unit definition: λ DNA (328 sites).

Reaction conditions:

66 mM potassium acetate, 33 mM Tris-acetate (pH 7.9), 10 mM magnesium acetate, 0.5 mM dithiothreitol, 0.1% Triton X-100, 100 μg/ml BSA. Incubate at **55°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 *Bsi*S I, >95% of the DNA fragments can be ligated and recut with this enzyme.

Heat inactivation: No.

Note: Blocked by overlapping *dcm* methylation.