

BshF I (Hae III)



Source: *Bacillus sphaericus*.

Cat.-No.	Size	Conc.
EN-110S	7,000 units	10 u/μl
EN-110L	35,000 units	10 u/μl

NB:
BSA is already included in the Buffer solution!

Buffer supplied: 10x B5 (incl. 10x BSA).

Substrate for unit definition: λ DNA (149 sites).

Reaction conditions:

50 mM potassium acetate, 20 mM Tris-acetate (pH 7.9), 10 mM magnesium acetate, 1 mM dithiothreitol, 100 μg/ml BSA.
Incubate at **37°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 *BshF I*, >95% of the DNA fragments can be ligated and recut with this enzyme.

Heat inactivation: 80°C for 20 minutes.

Note: Blocked by overlapping *dcm* methylation.