

# BamH I



Source: *Bacillus amyloliquefaciens* H.

---

Cat.-No.	Size	Conc.
EN-103S	7,500 units	10 u/μl
EN-103L	37,500 units	10 u/μl

**NB:**  
**BSA is already included in the Buffer solution!**

**Buffer supplied:** 10x *BamH I* (incl. 10x BSA).

**Substrate for unit definition:** λ DNA (5 sites)

**Reaction conditions:**

100 mM NaCl, 10 mM Tris-HCl (pH 7.9), 5 mM MgCl<sub>2</sub>, 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 50-fold overdigestion with *BamH I*, >95% of the DNA fragments can be ligated and recut with this enzyme.

**Star activity:**

Conditions of low ionic strength, high enzyme concentration, glycerol concentration >5%, or pH >8.0 may result in star activity.

**Heat inactivation:** 80°C for 20 minutes.