

# *PspP* I (*Sau96* I)



Source: *Psychrobacter immobilis* TA 137.

---

Cat.-No.	Size	Conc.
EN-126S	900 units	10 u/μl
EN-126L	4,500 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 (74 sites).

**Reaction conditions:**

50 mM NaCl, 10 mM Tris-HCl (pH 7.9), 10 mM MgCl<sub>2</sub>,  
1 mM dithiothreitol, 100 μg/ml BSA.  
Incubate at **25°C**.

**Storage buffer:**

50 mM NaCl, 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 *PspP* I, >95% of the  
DNA fragments can be ligated and recut with this  
enzyme.

**Note:** Incubation at 37°C results in 60% activity.

**Heat inactivation:** 55°C for 15 minutes.

Note: Blocked by overlapping *dcm* methylation.