



Source: *Nocardia rubra*.

Cat.-No.	Size	Conc.
EN-125S	700 units	10 u/μl
EN-125L	3,500 units	10 u/μl

Buffer supplied: 10x *Nru* I and 10x BSA.

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

Reaction conditions:

100 mM KCl, 50 mM Tris-HCl (pH 8.0), 10 mM MgCl₂, 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 *Nru* I, <20% of the DNA fragments can be ligated.

Star activity:

Large excess of the enzyme results in the appearance of star activity.

Heat inactivation: 65°C for 20 minutes.

Note: Blocked by overlapping *dam* methylation.