



Ava I

Cat. No. 25220-013
Conc.: 2-8 U/μl

Size: 300 units
Store at -20°C (not frost-free).



Description:

Ava I is isolated from a strain of *E. coli* that bears the cloned *Ava I* gene from *Anabaena variabilis*. One unit is the amount of enzyme required to cleave 1 μg of λ DNA in one hour at 37°C in the appropriate buffer.

Components:

25220-013 *Ava I*
Y92500 REACT[®] 2

Storage Buffer:

50 mM Tris-HCl (pH 7.5)
0.1 mM EDTA
10 mM 2-mercaptoethanol
500 μg/ml BSA
50% (v/v) glycerol
0.1% (w/v) Triton[®] X-100

Reaction Conditions:

Use REACT 2. REACT 2 is the 10X concentrate assay buffer supplied for use with *Ava I*. Dilute 1 part in 10 in final reaction mixture.

Final concentration:

50 mM Tris-HCl (pH 8.0)
10 mM MgCl₂
50 mM NaCl

Assay at 37°C.

Store buffer at 4°C or -20°C.

Doc. Rev.: 04/23/01

This product is distributed for laboratory research use only. CAUTION: Not for diagnostic use. The safety and efficacy of this product in diagnostic or other clinical uses has not been established.

For technical questions about this product, call the Life Technologies TECH-LINE[™] (800) 828-6686.

Quality Control:

Endonuclease Assay:

Cleavage only at recognition site at 100-fold excess digest of λ DNA.

Exonuclease Assay:

Using 1 pmol radiolabeled termini in a 50- μ l reaction and up to 20 units of enzyme for one hour, the following results were observed:

1 unit of enzyme removed \leq 0.3% label from 5' ends.

1 unit of enzyme removed \leq 0.3% label from 3' ends.

Ligation/Recut Assay:

On λ DNA: \geq 95% ligation in 16 hours at 22°C. 100% recut.

The enclosed buffers were assayed with the enzyme and met quality control specifications.

Note:

Partially resistant to heat inactivation (10 min, 65°C).

REFER TO THE GIBCO BRL CATALOGUE AND REFERENCE GUIDE FOR NOTES ON CONDITIONS WHICH AFFECT ENZYME ACTIVITY.

Triton[®] is a registered trademark of Rohm & Haas Co.