

# AluI

## Restriction enzyme for molecular biology applications

AluI is a restriction enzyme that recognizes AG<sup>CT</sup> sites.

10X Cutting Buffer is included that contains BSA, which enhances enzyme stability and binds to contaminants in DNA preps.

### Ordering Information

[Order Online »](#)

ENZ-GEN108-1000	1000U
-----------------	-------

### Manuals, SDS & CofA

[View Online »](#)

- High quality enzyme with stringent QC testing
- Includes 10X Cutting Buffer
- BSA premixed into buffers
- Suitable for molecular cloning, restriction site mapping, genotyping, Southern blotting, SNP



## Handling & Storage

Long Term Storage      -20°C

Shipping      Blue Ice

**Regulatory Status**      RUO - Research Use Only

## Product Details

**Activity**      One unit is defined as the amount of enzyme required to digest 1 µg of λ DNA in 1 hour at 37°C in a total reaction volume of 50 µl.

**Application Notes**      For molecular cloning, restriction site mapping, genotyping, Southern blotting, SNP.

**Concentration**      Alul 10,000 units/ml Includes 10X Cutting Buffer

**Formulation**      Liquid. In 100mM KCl, containing 10mM Tris-HCl, pH 7.4, 1mM DTT, 0.1mM EDTA, 50% glycerol and 200 µg/ml BSA.

After reconstitution, 1X Cutting Buffer: 50mM potassium acetate, 20mM Tris-acetate, 10mM magnesium acetate, 100 µg/ml BSA, pH 7.9, at 25°C.

**Quality Control**      **Exonuclease Activity (Radioactivity Release):**  
A 50 µl reaction in 1X Cutting Buffer containing 1 µg of a mixture of single and double stranded [<sup>3</sup>H] *E. coli* DNA and a minimum of 30 units of Alul incubated for 4 hours at 37°C releases <0.1% of the total radioactivity.

**Ligation and Recutting (Terminal Integrity):**  
After a 10-fold over digestion of λ DNA with Alul, ~75% of the DNA fragments can be ligated with T4 DNA ligase in 16 hours at 16°C. Of these ligated fragments, >95% can be recut with Alul.

**Non-Specific DNase Activity (16 Hour):**  
A 50 µl reaction in 1X Cutting Buffer containing 1 µg of λ DNA and a minimum of 50 Units of Alul incubated for 16 hours at 37°C results in a DNA pattern free of detectable nuclease degradation as determined by agarose gel electrophoresis.

**Reconstitution**      Reconstitute 10X Cutting Buffer with nuclease-free water.

**Source**      Produced in an *E. coli* strain that carries the cloned Alul gene from *Arthrobacter luteus*.



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)