

AluI

Restriction enzyme for molecular biology applications

AluI is a restriction enzyme that recognizes AG^{CT} 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 1µg of λ DNA in 1 hour at 37°C in a total reaction volume of 50µl.

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

Concentration Alul 10,000 units/mlIncludes 10X Cutting Buffer

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

After reconstitution, 1X Cutting Buffer: 50mM potassium acetate, 20mM Tris-acetate, 10mM magnesium acetate, 100µg/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 [³H] *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 AluI gene from *Arthrobacter luteus*.

Last modified: May 29, 2024



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-
be@enzolifesciences.com](mailto:info-be@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)