

# 176种内切酶，使用统一缓冲液，酶切如此简单！

## FastDigest快速内切酶

ThermoFisher  
SCIENTIFIC

| 内切酶名称               | 识别序列<br>5'→ 3'           | 反应温度<br>°C | 酶切时间 (min), 使用1 μL FastDigest内切酶 |                     |                           |                      | 完全酶切DNA<br>末端所需碱基对<br>数目 | 热失活          | 最长孵育时间<br>(h), 无星号<br>活性 | 甲基化敏感性               | 反应次数      | 货号       |
|---------------------|--------------------------|------------|----------------------------------|---------------------|---------------------------|----------------------|--------------------------|--------------|--------------------------|----------------------|-----------|----------|
|                     |                          |            | Lambda<br>DNA<br>1 μg/20 μL      | 质粒DNA<br>1 μg/20 μL | PCR产物<br>~0.2 μg/30<br>μL | 基因组DNA<br>1 μg/10 μL |                          |              |                          |                      |           |          |
| FastDigest AarI     | TTAATAA                  | 37         | 5                                | 5                   | 5                         | 5                    | 5                        | 65°C, 5 min  | 16                       | No effect            | 20        | FD2064   |
| FastDigest AasI     | GACNNNNJANGTC            | 37         | 5                                | 5                   | 5                         | 5                    | 1                        | 80°C, 10 min | 1                        | DS                   | 25        | FD1724   |
| FastDigest AatII    | GACGTC                   | 37         | 15                               | 20                  | 15                        | 15                   | 5                        | 80°C, 5 min  | 16                       | DS                   | 50        | FD0994   |
| FastDigest Acc65I   | GAGTACC                  | 37         | 5                                | 5                   | 5                         | 5                    | 3                        | 65°C, 5 min  | 16                       | DS DS                | 100       | FD0904   |
| FastDigest Adel     | CACNNN↓GTG               | 37         | 5                                | 5                   | 5                         | 5                    | 2                        | 80°C, 5 min  | 1                        | DS                   | 50        | FD1234   |
| FastDigest AjuI*    | ↓T/T(12GAAN)↓TTGG(11/6)↓ | 37         | 5                                | 5                   | 10                        | 5                    | ND                       | 65°C, 5 min  | 16                       | No effect            | 20        | FD1954   |
| FastDigest AluI     | AGCT                     | 37         | 15                               | 15                  | 15                        | 15                   | 4                        | 65°C, 5 min  | 16                       | No effect            | 100       | FD0014   |
| FastDigest Alw21I   | GWGCG↓C                  | 37         | 5                                | 5                   | 5                         | 5                    | 1                        | 80°C, 20 min | 16                       | No effect            | 100       | FD0024   |
| FastDigest Alw26I   | GTCTC(1/5)↓              | 37         | 5                                | 5                   | 5                         | 5                    | 1                        | 65°C, 5 min  | 16                       | DS                   | 100       | FD0034   |
| FastDigest Alw44I   | GATGCAC                  | 37         | 5                                | 60                  | 10                        | 5                    | 4                        | 80°C, 5 min  | 16                       | DS                   | 200       | FD0044   |
| FastDigest ApaI     | GGGCG↓C                  | 37         | 5                                | 5                   | 20                        | 10                   | 2                        | 65°C, 5 min  | 16                       | DS DS                | 300       | FD1414   |
| FastDigest BamHI    | G↓GATCC                  | 37         | 5                                | 5                   | 5                         | 5                    | 2                        | 80°C, 5 min  | 1                        | No effect            | 800/2,500 | FD0054/5 |
| FastDigest BclI     | T↓GATCA                  | 37         | 5                                | 5                   | 15                        | 5                    | 3                        | 80°C, 20 min | 16                       | DS                   | 300       | FD0724   |
| FastDigest BcnI     | CC↓ASGG                  | 37         | 5                                | 5                   | 5                         | 5                    | 3                        | 80°C, 20 min | 16                       | DS                   | 200       | FD0064   |
| FastDigest BclII    | A↓CTAGT                  | 37         | No sites                         | 5                   | 5                         | 5                    | 1                        | No           | 16                       | No effect            | 50        | FD1254   |
| FastDigest BclIII   | C↓ATRYAG                 | 37         | 5                                | 5                   | 5                         | 5                    | 2                        | 65°C, 20 min | 16                       | No effect            | 20        | FD1164   |
| FastDigest BfoI     | RGCGC↓Y                  | 37         | 5                                | 5                   | 5                         | 5                    | 2                        | 65°C, 10 min | 1                        | DS                   | 200       | FD2184   |
| FastDigest BglI     | GCCNNNNJANGGC            | 37         | 5                                | 5                   | 5                         | 5                    | 3                        | 65°C, 5 min  | 2                        | DS                   | 200       | FD0074   |
| FastDigest BglII    | A↓GATCT                  | 37         | 5                                | 20                  | 30                        | 20                   | 3                        | No           | 16                       | No effect            | 200       | FD0084   |
| FastDigest BmeI390I | CC↓ANGG                  | 37         | 5                                | 5                   | 5                         | 5                    | 4                        | No           | 16                       | DS DS                | 100       | FD1424   |
| FastDigest BmsI     | GCATC(5/9)↓              | 37         | 5                                | 5                   | 5                         | 5                    | 1                        | 65°C, 5 min  | 16                       | DS                   | 20        | FD2124   |
| FastDigest BoxI     | GACNNJANGTC              | 37         | 5                                | 5                   | 5                         | 5                    | 3                        | 80°C, 20 min | 16                       | DS                   | 100       | FD1434   |
| FastDigest BpiI     | G↓AGAC(2/6)↓             | 37         | 5                                | 5                   | 5                         | 5                    | 1                        | 65°C, 10 min | 16                       | No effect            | 20        | FD1014   |
| FastDigest BplI*    | ↓J(8/13GAG)N↓CTG(13/8)↓  | 37         | 15                               | 20                  | 30                        | 30                   | ND                       | 65°C, 5 min  | 16                       | No effect            | 20        | FD1314   |
| FastDigest Bpu10I   | CCTNAGC(↓5-2)↓           | 37         | 15                               | 15                  | 30                        | 15                   | 3                        | 80°C, 5 min  | 1                        | No effect            | 20        | FD1184   |
| FastDigest Bpu1102I | GC↓TNAGG                 | 37         | 5                                | 5                   | 5                         | 5                    | 2                        | 80°C, 5 min  | 16                       | No effect            | 50        | FD0094   |
| FastDigest BseDI    | C↓JGNGG                  | 37         | 5                                | 5                   | 5                         | 5                    | 3                        | 80°C, 5 min  | 16                       | No effect            | 20        | FD1084   |
| FastDigest BseGI    | GGATG(2/9)↓              | 37         | 5                                | 5                   | 5                         | 5                    | 2                        | 80°C, 5 min  | 16                       | No effect            | 100       | FD0874   |
| FastDigest BseII    | GATNNJANGTC              | 65         | 5                                | 5                   | 5                         | 5                    | 3                        | No           | 1                        | DS                   | 200       | FD1714   |
| FastDigest BseLI    | CCNNNNN↓ANGG             | 37         | 5                                | 5                   | 5                         | 5                    | 2                        | No           | 16                       | DS DS                | 100       | FD1204   |
| FastDigest BseMI    | GCAATG(2/9)↓             | 55         | 15                               | 15                  | 15                        | 15                   | 3                        | 80°C, 5 min  | 16                       | No effect            | 20        | FD1264   |
| FastDigest BseMII   | CTCAG(10/8)↓             | 55         | 15                               | 30                  | 15                        | 15                   | 2                        | 80°C, 5 min  | 16                       | No effect            | 20        | FD1404   |
| FastDigest BseNI    | ACTGG(1/-1)↓             | 65         | 5                                | 5                   | 5                         | 5                    | 2                        | 80°C, 5 min  | 16                       | No effect            | 100       | FD0884   |
| FastDigest BseSI    | GGGCG↓C                  | 37         | 5                                | 5                   | 5                         | 30                   | 3                        | 80°C, 15 min | 16                       | No effect            | 50        | FD1444   |
| FastDigest BseXI    | GCAGC(8/12)↓             | 65         | 15                               | 15                  | 15                        | 15                   | 3                        | 80°C, 15 min | 16                       | No effect            | 20        | FD1454   |
| FastDigest Bsh1236I | CG↓CG                    | 37         | 5                                | 5                   | 5                         | 5                    | 1                        | 80°C, 10 min | 16                       | DS                   | 100       | FD0924   |
| FastDigest Bsh1285I | CBRY↓CG                  | 37         | 15                               | 15                  | 15                        | 15                   | 3                        | 80°C, 15 min | 1                        | DS DS                | 100       | FD0894   |
| FastDigest BshNI    | G↓GYRCC                  | 37         | 5                                | 60                  | 5                         | 15                   | 2                        | 65°C, 10 min | 16                       | DS DS                | 300       | FD1004   |
| FastDigest BshTl    | A↓CCGGT                  | 37         | 5                                | 5                   | 5                         | 5                    | 3                        | 80°C, 5 min  | 16                       | DS DS                | 20        | FD1464   |
| FastDigest Bsp119I  | TT↓GCGAA                 | 37         | 5                                | 5                   | 5                         | 5                    | 1                        | 80°C, 5 min  | 16                       | DS                   | 200       | FD0124   |
| FastDigest Bsp120I  | G↓GGCCC                  | 37         | 5                                | 5                   | 15                        | 5                    | 3                        | 80°C, 10 min | 16                       | DS DS                | 200       | FD0134   |
| FastDigest Bsp1407I | T↓GTACA                  | 37         | 5                                | 5                   | 5                         | 5                    | 3                        | 80°C, 10 min | 16                       | No effect            | 100       | FD0934   |
| FastDigest Bsp143I  | J↓GATC                   | 37         | 5                                | 5                   | 10                        | 5                    | 4                        | 65°C, 20 min | 16                       | DS                   | 40        | FD0784   |
| FastDigest BspLI    | GGN↓AAC                  | 37         | 5                                | 5                   | 5                         | 5                    | 2                        | 65°C, 20 min | 16                       | DS DS                | 20        | FD1154   |
| FastDigest BspOI    | GCTAG↓C                  | 37         | 5                                | 5                   | 15                        | 5                    | 3                        | 80°C, 5 min  | 1                        | No effect            | 20        | FD2044   |
| FastDigest BspTl    | C↓TTAAG                  | 37         | 5                                | 5                   | 5                         | 5                    | 4                        | No           | 16                       | No effect            | 150       | FD0834   |
| FastDigest Bst1107I | GTA↓TAC                  | 37         | 5                                | 5                   | 5                         | 10                   | 3                        | No           | 6                        | DS                   | 100       | FD0704   |
| FastDigest BstXI    | CCANNNNNJANGTG           | 37         | 5                                | 5                   | 5                         | 5                    | 4                        | 80°C, 5 min  | 0.5                      | DS DS                | 100       | FD1024   |
| FastDigest Bsu15I   | AT↓CGAT                  | 37         | 5                                | 5                   | 5                         | 5                    | 2                        | 65°C, 15 min | 16                       | DS DS                | 100       | FD0144   |
| FastDigest BsuRI    | GG↓CC                    | 37         | 5                                | 5                   | 5                         | 5                    | 3                        | No           | 16                       | No effect            | 400       | FD0154   |
| FastDigest BveI     | ACCTGC(4/8)↓             | 37         | 15                               | 15                  | 60                        | 20                   | 5                        | 65°C, 5 min  | 16                       | DS                   | 50        | FD1744   |
| FastDigest CaiI     | CAGNNN↓CTG               | 37         | 5                                | 5                   | 5                         | 5                    | 4                        | 65°C, 10 min | 16                       | DS DS                | 50        | FD1394   |
| FastDigest Cfr10I   | R↓CCGGY                  | 37         | 5                                | 5                   | 10                        | 5                    | 2                        | No           | 2                        | DS                   | 50        | FD0184   |
| FastDigest Cfr13I   | G↓GNCC                   | 37         | 5                                | 5                   | 5                         | 5                    | 2                        | No           | 16                       | DS DS                | 100       | FD0194   |
| FastDigest CpOI     | CG↓GGWCCG                | 37         | 5                                | 5                   | 5                         | 5                    | 1                        | 65°C, 5 min  | 16                       | DS                   | 20        | FD0744   |
| FastDigest CseI     | GACGG(5/10)↓             | 37         | 15                               | 15                  | 30                        | 15                   | 2                        | 80°C, 5 min  | 16                       | DS                   | 20        | FD1904   |
| FastDigest CsiI     | A↓CCWGGT                 | 37         | 5                                | 10                  | 5                         | 15                   | 4                        | 65°C, 5 min  | 16                       | No effect            | 50        | FD2114   |
| FastDigest Csp6I    | G↓TAC                    | 37         | 5                                | 5                   | 5                         | 5                    | 2                        | 80°C, 10 min | 16                       | No effect            | 100       | FD0214   |
| FastDigest DpnI     | Gm↓AATC                  | 37         | ND                               | 5                   | Not applicable            | 10                   | 1                        | 80°C, 5 min  | 16                       | Does not cut dam DNA | 100       | FD1704   |
| FastDigest DraI     | TTT↓AAA                  | 37         | 5                                | 5                   | 5                         | 5                    | 4                        | 65°C, 5 min  | 16                       | No effect            | 200       | FD0224   |
| FastDigest Eam1104I | CTCTCT(1/4)↓             | 37         | 5                                | 5                   | 5                         | 5                    | 3                        | 80°C, 5 min  | 6                        | No effect            | 50        | FD0234   |
| FastDigest Eam1105I | GACNNN↓ANGTC             | 37         | 5                                | 5                   | 5                         | 5                    | 3                        | 65°C, 5 min  | 1                        | No effect            | 100       | FD0244   |
| FastDigest Ecl136II | GAG↓CTC                  | 37         | 5                                | 5                   | 5                         | 5                    | 1                        | 65°C, 5 min  | 6                        | DS                   | 100       | FD0254   |
| FastDigest Eco105I  | TAC↓GTA                  | 37         | 5                                | 5                   | 5                         | 5                    | 3                        | 65°C, 5 min  | 16                       | DS                   | 50        | FD0404   |
| FastDigest Eco130I  | C↓CWWGG                  | 37         | 5                                | 5                   | 20                        | 20                   | 3                        | 65°C, 5 min  | 16                       | No effect            | 200       | FD0414   |
| FastDigest Eco147I  | AGG↓CCT                  | 37         | 5                                | 5                   | 5                         | 5                    | 3                        | 80°C, 10 min | 16                       | DS                   | 100       | FD0424   |
| FastDigest Eco31I   | GGTCTC(1/5)↓             | 37         | 5                                | 5                   | 5                         | 5                    | 3                        | 65°C, 5 min  | 16                       | DS DS                | 100       | FD0294   |
| FastDigest Eco32I   | GAT↓ATC                  | 37         | 5                                | 5                   | 5                         | 5                    | 2                        | No           | 16                       | No effect            | 400       | FD0304   |
| FastDigest Eco47I   | G↓GWCC                   | 37         | 5                                | 5                   | 5                         | 5                    | 1                        | 80°C, 20 min | 16                       | DS DS                | 200       | FD0314   |
| FastDigest Eco47III | AGC↓GCT                  | 37         | 5                                | 5                   | 5                         | 5                    | 4                        | 65°C, 5 min  | 16                       | DS                   | 20        | FD0324   |
| FastDigest Eco52I   | C↓GGCCG                  | 37         | 5                                | 20                  | 20                        | 5                    | 3                        | 65°C, 5 min  | 16                       | DS                   | 50        | FD0334   |
| FastDigest Eco57I   | CTGAAG(16/14)↓           | 37         | 15                               | 30                  | 15                        | 15                   | 2                        | 65°C, 5 min  | 16                       | No effect            | 20        | FD0344   |
| FastDigest Eco72I   | CAC↓GTG                  | 37         | 5                                | 5                   | 5                         | 5                    | 2                        | 80°C, 10 min | 1                        | DS                   | 200       | FD0364   |
| FastDigest Eco81I   | CC↓TNAGG                 | 37         | 5                                | 5                   | 5                         | 5                    | 2                        | 80°C, 10 min | 16                       | No effect            | 50        | FD0374   |
| FastDigest Eco88I   | C↓YCGRG                  | 37         | 5                                | 5                   | 5                         | 5                    | 2                        | 65°C, 5 min  | 16                       | DS                   | 200       | FD0384   |
| FastDigest Eco91I   | G↓GTNACC                 | 37         | 5                                | 5                   | 5                         | 5                    | 2                        | 65°C, 10 min | 2                        | No effect            | 200       | FD0394   |
| FastDigest EcoN109I | RG↓GNCYC                 | 37         | 5                                | 10                  | 15                        | 5                    | 2                        | 65°C, 5 min  | 16                       | DS                   | 200       | FD0264   |
| FastDigest EcoRI    | G↓AATTC                  | 37         | 5                                | 5                   | 20                        | 5                    | 2                        | 80°C, 5min   | 0.5                      | DS DS                | 800/2,500 | FD0274/5 |
| FastDigest EheI     | GGC↓GGC                  | 37         | 5                                | 5                   | 5                         | 5                    | 2                        | 65°C, 5 min  | 6                        | DS                   | 50        | FD0444   |
| FastDigest Esp3I    | CGTCTC(1/5)↓             | 37         | 5                                | 5                   | 5                         | 5                    | 1                        | 65°C, 10 min | 6                        | DS                   | 20        | FD0454   |
| FastDigest FagI     | GGGAC(10/14)↓            | 37         | 15                               | 15                  | 15                        | 15                   | 3                        | 65°C, 5 min  | 16                       | DS                   | 20        | FD1814   |
| FastDigest FokI     | GGATG(9/13)↓             | 37         | 5                                | 5                   | 5                         | 5                    | 1                        | 65°C, 5 min  | 1                        | No effect            | 100       | FD2144   |
| FastDigest FspAI    | RTGC↓GAY                 | 37         | 5                                | 5                   | 5                         | 5                    | 4                        | 65°C, 5 min  | 16                       | DS                   | 20        | FD1664   |
| FastDigest FspBI    | C↓TAG                    | 37         | 5                                | 151                 |                           |                      |                          |              |                          |                      |           |          |