

pCBG99-Basic Vector Restriction Enzyme Tables

The following restriction enzyme tables were constructed using DNASTAR® sequence analysis software. This information has not been verified by restriction digestion with each enzyme listed. The location given specifies the 3'-end of the cut DNA (the base to the left of the cut site). Vector sequences are also available in the GenBank® database (GenBank®/EMBL Accession Number AY258595) and on the Internet at: www.promega.com/vectors/

Restriction Enzymes That Cut the pCBG99-Basic Vector 1-5 Times.

| Enzyme | # of Sites | Location | Enzyme | # of Sites | Location |
|-----------------|------------|------------------------------|----------------|------------|------------------------------|
| AatII | 1 | 314 | Eco52I | 3 | 1730, 1734, 4626 |
| AccB7I | 1 | 138 | EcoICRI | 1 | 9 |
| AccI | 2 | 794, 1986 | FseI | 1 | 1736 |
| Acc65I | 1 | 1 | FspI | 2 | 3350, 4523 |
| AcyI | 2 | 311, 3665 | HaeII | 5 | 1121, 2113, 2483, 4122, 4130 |
| AflIII | 4 | 15, 753, 1590, 2235 | HgaI | 5 | 718, 2346, 2924, 3654, 4055 |
| Alw26I | 3 | 859, 3189, 3965 | HincII | 4 | 399, 1635, 1877, 1987 |
| Alw44I | 2 | 2549, 3795 | HindIII | 4 | 399, 1635, 1877, 1987 |
| AlwNI | 1 | 2651 | HindIII | 1 | 53 |
| ApaI | 1 | 1228 | HpaI | 1 | 1877 |
| AspHI | 5 | 11, 177, 2553, 3714, 3799 | Hsp92I | 2 | 311, 3665 |
| AvaI | 3 | 26, 32, 1135 | KpnI | 1 | 5 |
| AvaII | 2 | 3266, 3488 | MluI | 2 | 15, 753 |
| BamHI | 1 | 1979 | NciI | 5 | 27, 28, 2615, 3311, 3662 |
| BanII | 5 | 11, 33, 973, 1228, 4206 | NcoI | 1 | 86 |
| BbsI | 3 | 453, 989, 2064 | NheI | 1 | 21 |
| BclI | 2 | 882, 1400 | NotI | 1 | 4626 |
| BglI | 2 | 3248, 4516 | NruI | 2 | 1205, 1472 |
| BglII | 1 | 36 | NsiI | 1 | 1468 |
| BsaI | 2 | 859, 3189 | NspI | 1 | 2239 |
| BsaAI | 1 | 4277 | PaeR7I | 1 | 32 |
| BsaBI | 1 | 1978 | PflMI | 1 | 138 |
| BsaHI | 2 | 311, 3665 | Ppu10I | 1 | 1464 |
| Bsp120I | 1 | 1224 | PshAI | 1 | 2050 |
| BspHI | 2 | 2955, 3963 | PspAI | 1 | 26 |
| BspMI | 1 | 4756 | PvuI | 2 | 3498, 4544 |
| BsrBRI | 1 | 1978 | PvuII | 2 | 768, 1432 |
| BssSI | 3 | 976, 2408, 3792 | SacI | 1 | 11 |
| BstXI | 1 | 292 | Sall | 1 | 1985 |
| BstZI | 3 | 1730, 1734, 4626 | Scal | 3 | 1609, 3608, 4691 |
| ClaI | 3 | 1972, 4684, 4788 | SinI | 2 | 3266, 3488 |
| DraI | 5 | 621, 1938, 2994, 3013, 3705 | SmaI | 1 | 28 |
| DraII | 1 | 1258 | SrfI | 1 | 28 |
| DraIII | 2 | 1006, 4280 | SspI | 5 | 477, 555, 3932, 4485, 4600 |
| DrdI | 2 | 2343, 4324 | StyI | 3 | 86, 206, 1301 |
| DsaI | 1 | 86 | VspI | 1 | 3300 |
| EaeI | 5 | 1055, 1730, 1734, 3516, 4626 | XbaI | 1 | 1717 |
| EagI | 3 | 1730, 1734, 4626 | XhoI | 1 | 32 |
| EarI | 3 | 2119, 3923, 4561 | XmaI | 1 | 26 |
| EclHKI | 1 | 3128 | XmnI | 1 | 3727 |
| Eco47III | 2 | 1119, 2111 | | | |

Restriction Enzymes That Do Not Cut the pCBG99-Basic Vector.

| | | | | | | | |
|---------------|----------------|---------------|---------------|-------------|--------------|----------------|------|
| AccIII | BbrPI | Bst98I | EcoNI | NdeI | PstI | SpeI | XcmI |
| AflII | BbuI | BstEII | EcoRI | Pacl | RsrII | SphI | |
| AgeI | BlpI | Bsu36I | EcoRV | PinAI | SacII | SpII | |
| AscI | Bpu1102I | CspI | EheI | PmeI | SfiI | Sse8387I | |
| AvrII | BsrGI | Csp45I | I-PpoI | PmlI | SgfI | StuI | |
| Ball | BssHIII | Eco72I | KasI | PpuMI | SgrAI | Swal | |
| BbeI | Bst1107I | Eco81I | NarI | Psp5II | SnaBI | Tth111I | |

Note: The enzymes in boldface type are available from Promega.