## **Digestion of agarose-embedded DNA**

Pub. No. MAN0011974 Rev. B.0

Restriction digestion of DNA embedded in agarose is a technique used for analysis of extremely large DNA fragments. Complete digestion of substrate DNA is affected by substrate purity, units of enzyme, and incubation time. The following protocol describes a general procedure for the digestion of 1 µg of agarose-embedded DNA in 16 hours.

## Digest agarose-embedded DNA with restriction enzyme

- 1. Embed the substrate DNA in 1% low melting temperature agarose at a concentration of 1 µg DNA / 30 µL agarose.
- 2. Prepare ~30 µL agarose plugs with a gel syringe dispenser or similar agarose dispensing system.
- 3. Equilibrate the plug in 100  $\mu$ L of the appropriate 1X restriction enzyme buffer for 15 minutes.
- 4. Place the plug in 100 µL of fresh 1X buffer containing the restriction enzyme (see Table 1 for recommended number of units).
- 5. Incubate at 37°C for 16 hours.

Note: Certain enzymes may require incubation at higher or lower temperatures. Refer to Table 1 for details on specific enzymes.

## Table 1. Recommended units of enzyme for digestion of agarose-embeded DNA

Enzyme	Cat. no .	Units of enzyme required for digestion in 16 hours <sup>[1]</sup>
Aanl (Psil)	ER2061	5
Aarl	ER1581	5
AatII	ER0991	5
Acc65I (Asp718I)	ER0901	5
Adel (Dralli)	ER1231	10
Ajil (BmgBl)	ER1941	5
Ajul	ER1951	5
Alw44I	ER0041	5
Apal	ER1411	5
BamHI	ER0051	5
Baul (BssSI)	ER1841	5
Bcul (Spel)	ER1251	10
Bful (BciVI)	ER1501	5
Bgll	ER0071	5
BgllI	ER0081	5
Boxl (PshAl)	ER1431	10
Bpil (Bbsl)	ER1011	5
Bpll	ER1311	10



Enzyme	Cat. no .	Units of enzyme required for digestion in 16 hours <sup>[1]</sup>
Bpu1102I (BlpI)	ER0091	5
BseSI (Bme1580I)	ER1441	5
Bsh1285I (BsiEI)	ER0891	5
BshNI (BanI)	ER1001	5
BshTl (Agel)	ER1461	20
Bsp68I (Nrul)	ER0111	5
Bsp119I (BstBI)	ER0121	5
Bsp120I (Psp0MI)	ER0131	5
Bsp1407I (BsrGI)	ER0931	10
Bsp0I (BmtI)	ER2041	20
BspTI (AflII)	ER0831	5
Bst1107I (BstZ17I)	ER0701	5
BstXI	ER1021	5
Bsu15I (ClaI)	ER0141	5
Cail (AlwNI)	ER1391	5
Cfr9l (Xmal)	ER0171	5
Cfr10l (BsrFl)	ER0181	5
Cfr42I (SacII)	ER0201	20
Dral	ER0221	5
Eam1104I (Earl)	ER0231	5
Eam1105I (AhdI)	ER0241	5
Ecl136II (EcoICRI)	ER0251	5
Eco24I (BanII)	ER0281	5
Eco31I (Bsal)	ER0291	5
Eco32I (EcoRV)	ER0301	5
Eco47III (Afel)	ER0321	5
Eco52I (EagI)	ER0331	5
Eco72I (PmlI)	ER0361	5
Eco811 (Bsu361)	ER0371	5
Eco88I (Aval)	ER0381	5
Eco91I (BstEII)	ER0391	5
Eco105I (SnaBI)	ER0401	5
Eco130I (Styl)	ER0411	5
Eco147I (Stul)	ER0421	5
EcoO109I (Drall)	ER0261	5
EcoRI	ER0271	5
Ehel (Sfol)	ER0441	20

Enzyme	Cat. no .	Units of enzyme required for digestion in 16 hours <sup>[1]</sup>
Esp3I (BsmBI)	ER0451	5
FspBI (Bfal)	ER1761	10
Hin1I (BsaHI)	ER0471	5
HindIII	ER0501	5
Kpnl	ER0521	5
Kpn2l (BspEl)	ER0531	5 (55°C)
KspAl (Hpal)	ER1031	5
Lgul (Sapl)	ER1931	5
MauBI	ER2081	5
Mbil (BsrBl)	ER1271	5
MlsI (MscI)	ER1211	20
Mlul	ER0561	5
Mssl (Pmel)	ER1341	5
Munl (Mfel)	ER0751	5
Mva1269I (Bsml)	ER0961	5
Ncol	ER0571	5
Ndel	ER0581	5
Nhel	ER0971	5
Notl	ER0591	5
Nsbl (Fspl)	ER1221	5
Olil (Alel)	ER1631	5
Pael (Sphl)	ER0601	5
Pacl	ER2201	5
Pagl (BspHI)	ER1281	5
Pasl	ER1861	10 (55°C)
Paul (BssHII)	ER1091	10
Pdil (Nael)	ER1521	20
Pdml (Xmnl)	ER1531	10
Pfl23II (BsiWI)	ER0851	5
Pfol	ER1751	10
Ppu21I (BsaAI)	ER1971	5
Pscl (Pcil)	ER1871	5
Psp5II (PpuMI)	ER0761	5
Psp1406I (AclI)	ER0941	10
Pstl	ER0611	5
Psyl (Tth111I)	ER1331	5
Pvul	ER0621	5

Enzyme	Cat. no .	Units of enzyme required for digestion in 16 hours <sup>[1]</sup>
Pvull	ER0631	5
Rsel (Msll)	ER2001	5
Sacl	ER1131	5
Sall	ER0641	5
Scal	ER0431	20
Sdal (Sbfl)	ER1191	5
SfaAl (AsiSI)	ER2091	5
Sfil	ER1821	5 (50°C)
Sgel	ER2211	3
SgrDl	ER2031	5
Sgsl (Ascl)	ER1891	5
Smal	ER0661	5 (30°C)
Smil (Swal)	ER1241	10 (30°C)
Smol (Smll)	ER1981	5 (55°C)
Sspl	ER0771	5
SspDI (Kasl)	ER2191	5
Van91I (PflMI)	ER0711	5
Vspl (Asel)	ER0911	5
Xagl (EcoNI)	ER1301	5
Xapl (Apol)	ER1381	5
Xbal	ER0681	5
Xhol	ER0691	5
XmaJI (AvrII)	ER1561	10
Xmil (Accl)	ER1481	5
I-Scel	ER1771	20

<sup>[1]</sup> Unless otherwise noted in parentheses, incubation temperature is 37°C.

The information in this guide is subject to change without notice.

## DISCLAIMER

TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

Corporate entity: Life Technologies Corporation | Carlsbad, CA 92008 USA | Toll Free in USA 1 800 955 6288

©2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

For support visit thermofisher.com/support or email techsupport@lifetech.com



thermofisher.com