

Find the NEB Exonuclease Best Suited for your Application



Potential Substrates

DOUBLE-STRANDED



Supercoiled



Nicked circular*



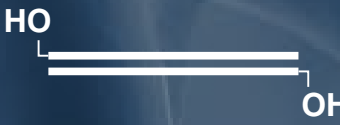
SINGLE-STRANDED



Circular



What do I want for ending material(s)**



Degraded in the presence of Mg²⁺



Complete degradation with long incubation



Preferentially degrades phosphorylated 5' strand



Recommended Exonuclease

Exonuclease T	M0265S/L
Exonuclease I (<i>E. coli</i>)	M0293S/L
Thermolabile Exonuclease I	M0568S/L
<i>Msz</i> Exonuclease I	M0527S
RecJ _f	M0264S/L

Exonuclease VII	M0379S/L
-----------------	----------

Nuclease P1 [†]	M0660S
--------------------------	--------

Exonuclease VIII, truncated	M0545S/L
-----------------------------	----------

Exonuclease V (RecBCD)	M0345S/L
------------------------	----------

Lambda Exonuclease	M0262S/L
--------------------	----------

Exonuclease III (<i>E. coli</i>) ^{††}	M0206S/L
--	----------

T7 Exonuclease	M0263S/L
----------------	----------

T5 Exonuclease	M0663S/L
----------------	----------

* Nicked plasmid is prepared by treating supercoiled DNA with a nicking enzyme.
** End products less than 10-mer can be removed by Monarch® cleanup kits and are not shown.
† Although a single-strand specific nuclease (ssDNA and RNA-specific), Nuclease P1 does display some activity toward dsDNA in Nuclease P1 Reaction Buffer. If preferentially degrading single-stranded nucleic acids (ssDNA or RNA) in the presence of double stranded DNA (dsDNA), we recommend using Nuclease P1 in 1X NEBuffer® r1.1 to limit activity on dsDNA while maintaining single-strand nuclease activity.
†† Extended incubation or excess enzyme may results in ssDNA deviation.