

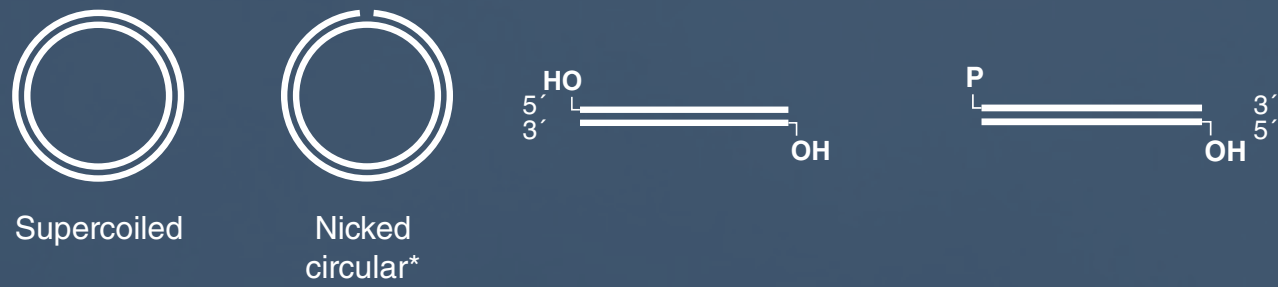
# Find the NEB Exonuclease Best Suited for your Application



be INSPIRED  
drive DISCOVER  
stay GENUINE

## DOUBLE-STRANDED

## SINGLE-STRANDED



### What do I want for ending material(s)?†

Starting Material	Exonuclease	NEB #
Supercoiled, Nicked circular, Linear (5' OH, 3' OH), Linear (5' P, 3' OH)	Exonuclease T	M0265S/L
Supercoiled, Nicked circular, Linear (5' OH, 3' OH), Linear (5' P, 3' OH)	Exonuclease I ( <i>E. coli</i> )	M0293S/L
Supercoiled, Nicked circular, Linear (5' OH, 3' OH), Linear (5' P, 3' OH)	Thermolabile Exonuclease I	M0568S/L
Supercoiled, Nicked circular, Linear (5' OH, 3' OH), Linear (5' P, 3' OH)	Exonuclease VII	M0379S/L
Supercoiled, Nicked circular, Linear (5' OH, 3' OH), Linear (5' P, 3' OH)	Degraded by magnesium	
Supercoiled, Nicked circular, Linear (5' OH, 3' OH), Linear (5' P, 3' OH)	Nuclease P1	M0660S
Supercoiled, Nicked circular, Linear (5' OH, 3' OH), Linear (5' P, 3' OH)	Exonuclease VIII, truncated	M0545S/L
Supercoiled, Nicked circular, Linear (5' OH, 3' OH), Linear (5' P, 3' OH)	Exonuclease V (RecBCD)	M0345S/L
Supercoiled, Nicked circular, Linear (5' OH, 3' OH), Linear (5' P, 3' OH)	Lambda Exonuclease	M0262S/L
Supercoiled, Nicked circular, Linear (5' OH, 3' OH), Linear (5' P, 3' OH)	Exonuclease III ( <i>E. coli</i> )	M0206S/L
Supercoiled, Nicked circular, Linear (5' OH, 3' OH), Linear (5' P, 3' OH)	T7 Exonuclease	M0263S/L
Supercoiled, Nicked circular, Linear (5' OH, 3' OH), Linear (5' P, 3' OH)	T5 Exonuclease	M0363S/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.