

Dephosphorylation of Proteins

This protocol is for the Dephosphorylation of Proteins

Reaction Mixture:

1X Thermo Scientific™ FastAP™ reaction buffer,
0.1 - 0.2 mg/mL of phosphoprotein,
10 u of FastAP Thermosensitive Alkaline Phosphatase.
Incubate at 37 °C for 1 hour.

For example: if you are doing a 20 µL reaction setup you need 2 µL 10x FastAP buffer, 2-4 µg of protein (to be in the range of 0.1 - 0.2 mg/mL) and 10 U of FastAP Thermosensitive Alkaline Phosphatase (1 u/µL).

Note:

- The reaction can be stopped by addition of a final concentration of 50 mM EDTA (#R1021) or by addition of a final concentration of 10 mM sodium orthovanadate (Na₃VO₄).
- The optimal incubation time and the enzyme concentration must be determined experimentally for each substrate

www.thermoscientific.com/onebio

© 2013 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

Europe

Customer Service
cs.molbio.eu@thermofisher.com

Technical Support
ts.molbio.eu@thermofisher.com

Tel 00800 222 00 888
Fax 00800 222 00 889

United States

Customer Service
cs.molbio@thermofisher.com

Technical Support
ts.molbio@thermofisher.com

Tel 800 235 9880
Fax 800 292 6088

Canada

Customer Service
cs.molbio@thermofisher.com

Technical Support
ts.molbio@thermofisher.com

Tel 800 340 9026
Fax 800 472 8322

Thermo
S C I E N T I F I C

Part of Thermo Fisher Scientific