

DnaK (E. coli) monoclonal antibody (8E2/2)

DnaK possesses an autophosphorylation activity and a weak 5-nucleotidase activity, cleaving the 5 phosphate groups from both ribose and deoxyribose nucleotides. The role of DnaK in ATP-dependent protein-protein interactions has been extended to normal *E. coli* physiology, where, like eukaryotic Hsp70 homologs, it is thought to participate in the assembly/disassembly of protein complexes.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 7

[View Online »](#)

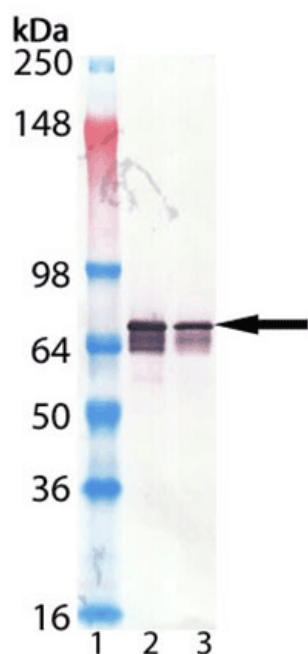
Ordering Information

[Order Online »](#)

ENZ-ABS680-0200	200µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western Blot analysis of ENZ-ABS680: Lane 1: MW Marker, Lane 2: DnaK (*E. coli*) recombinant protein (Prod. No. ADI-SPP-630), Lane 3: *E. coli* cell lysate.

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Alternative Name	Hsp70
Application	WB
Application Notes	Detects a band of ~70kDa by Western blot.
Clone	8E2/2
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.
Host	Mouse
Immunogen	E. coli DnaK.
Isotype	IgG1
Purity Detail	Protein G affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:10,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from hybridoma tissue culture supernatant.
Species Reactivity	E. coli
UniProt ID	P0A6Y8 (strain K12)
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)