

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: 71

[View Online »](#)

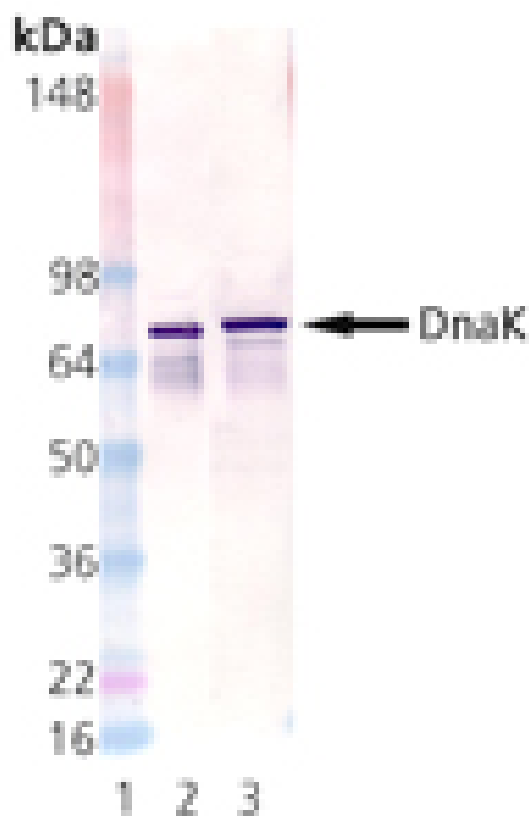
Ordering Information

[Order Online »](#)

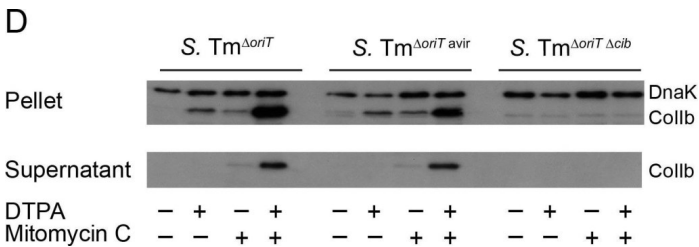
ADI-SPA-880-J	1mg
ADI-SPA-880-D	50µg
ADI-SPA-880-F	200µg

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis: Lane 1: MW Marker, Lane 2: DnaK active recombinant protein (Prod. No. ADI-SPP-630), Lane 3: E. coli cell lysate.



Expression of *S. Tm* Collb is induced by iron limitation and the SOS response. (A) Organization of the Collb locus showing location of the Fur and LexA repressor binding sites in the Collb (*cib*) promoter and the immunity protein gene (*imm*). Overnight cultures of *S. Tm*^{ΔoriT} (B) and *S. Tm*^{ΔoriT avir} (C) carrying the reporter plasmid *pcib-luc* were re-inoculated 1:20 in fresh LB with indicated supplements (0.25 μg/ml mitomycin C; 100 mM DTPA) and grown under aeration for 4 h. Cultures were normalized to OD600, bacteria were harvested and luciferase-activity was measured in bacterial lysates. Relative luminescence units (RLU) per unit OD600 are indicated. (D) Overnight cultures of indicated *S. Tm* strains were re-inoculated 1:20 in fresh LB with indicated supplements (0.25 μg/ml mitomycin C; 100 mM DTPA) and grown under aeration for 4 h. OD600 of the cultures was normalized, bacteria were harvested and Collb was detected in bacterial lysates as well as in the culture supernatant by immunoblot using an affinity-purified rabbit-α-Collb antiserum. *S. Tm* DnaK was detected as loading control.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Inflammation fuels colicin Ib-dependent competition of *Salmonella* serovar Typhimurium and *E. coli* in enterobacterial blooms. *PLoS Pathog* (2014)

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

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.

GenBank ID AP009048

Host Mouse

Immunogen E. coli DnaK.

Isotype IgG1

Purity Detail Protein G affinity purified.

Recommendation 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 ascites.

Species Reactivity E. coli

UniProt ID

P0A6Y8 (strain K12)



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee.