

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 »](#)

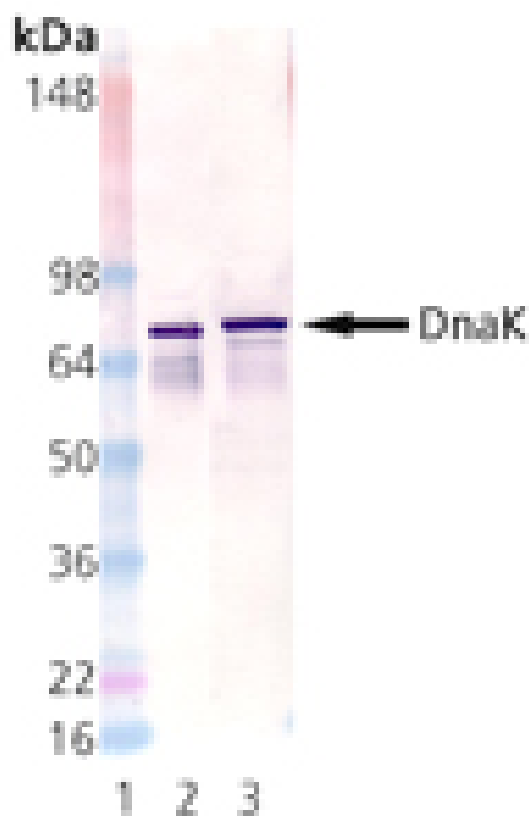
Ordering Information

[Order Online »](#)

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

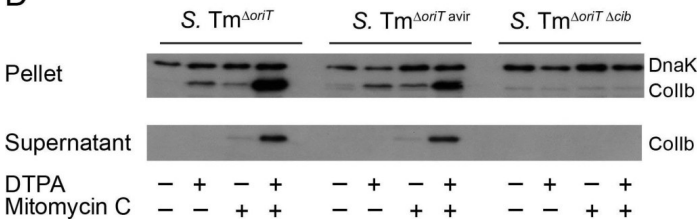
Manuals, SDS & CofA

[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.

D



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* Δ oriT (B) and *S. Tm* Δ oriT avir (C) carrying the reporter plasmid *pcib-luc* were re-inoculated 1:20 in fresh LB with indicated supplements (0.25 μ g/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 μ g/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 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 ascites.
Species Reactivity	E. coli
UniProt ID	P0A6Y8 (strain K12)

Worry-free Guarantee

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

Last modified: May 29, 2024



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)