

DnaK (E. coli), (recombinant)

Recombinant DnaK (E. coli) that functions as a molecular chaperone, assisting protein folding and complex assembly.

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.

Citations: 9

[View Online »](#)

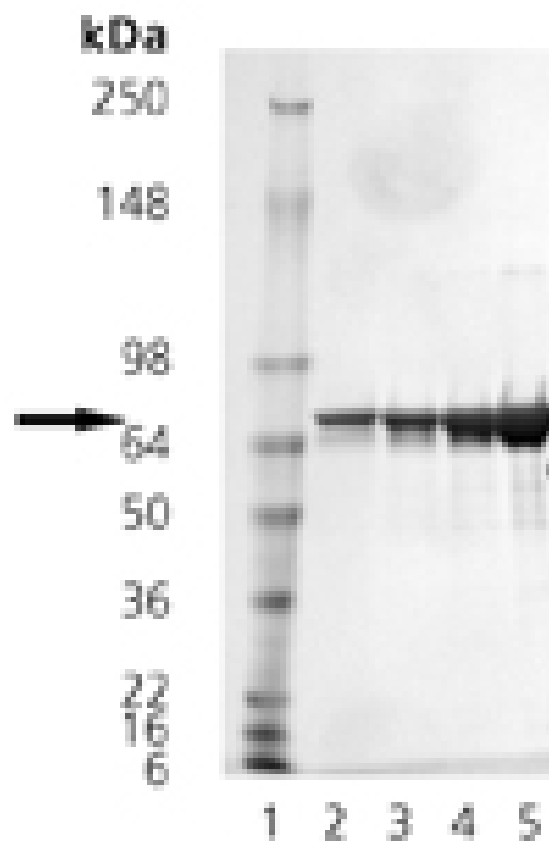
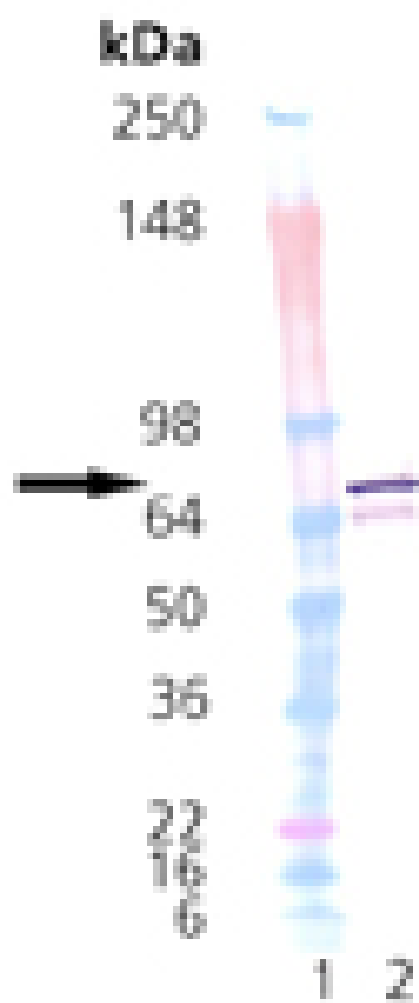
Ordering Information

[Order Online »](#)

ADI-SPP-630-J	1mg
ADI-SPP-630-D	50µg
ADI-SPP-630-F	200µg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name Heat shock protein 70, HSP70

Application Notes ATPase activity assay (positive). Western blot control.

Formulation Liquid. In 40mM TRIS, pH 7.5, containing 80mM sodium chloride, 0.8mM DTT, 0.08mM PMSF, and 20% glycerol.

MW ~70kDa

Purity ≥95% (SDS-PAGE; Western blot)

Purity Detail Purified by multi-step chromatography.

Source Produced in *E. coli*.

UniProt ID P0A6Y8 (strain K12)

Last modified: October 9, 2025



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)