

PRODUCT INFORMATION



Hsp40 (human recombinant)

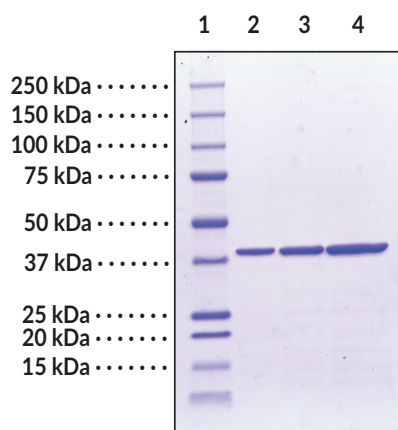
Item No. 23406

Overview and Properties

Synonym:	Heat Shock Protein 40
Source:	Recombinant, full-length, N-terminal histidine-tagged Hsp40 protein purified from <i>E. coli</i>
Amino acids:	2-340 (full length)
Uniprot No.:	P25685
Molecular Weight:	40.1 kDa
Storage:	-80°C (as supplied); avoid freeze/thaw cycles by storing protein in aliquots
Stability:	≥1 year
Purity:	<i>batch specific</i> (≥90% estimated by SDS-PAGE)
Supplied in:	50 mM HEPES, pH 8.0, 150 mM sodium chloride, and 10% glycerol
Protein	
Concentration:	<i>batch specific</i> mg/ml

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Image



Lane 1: MW Markers
Lane 2: Hsp40 (1 µg)
Lane 3: Hsp40 (2 µg)
Lane 4: Hsp40 (4 µg)

Representative gel image shown; actual purity may vary between each batch.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/09/2018

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Heat shock protein 40 (Hsp40) is an Hsp70 co-chaperone protein that regulates the ATP-dependent binding of Hsp70 (Item Nos. 23002 | 22739) to target proteins.¹ There are 44 members of the Hsp40 family in humans, and each Hsp40 contains a unique polypeptide-binding domain that binds and delivers specific target proteins to Hsp70. Hsp40 also contains a conserved HPD tripeptide that is critical to its role in driving conversion of Hsp70 from its ATP-bound form to its ADP-bound form to facilitate target loading.

Reference

1. Fan, C.-Y., Lee, S., and Cyr, D.M. Mechanisms for regulation of Hsp70 function by Hsp40. *Cell Stress Chaperones* **8(4)**, 309-316 (2003).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM