PRODUCT INFORMATION



Hsc70 (human, recombinant)

Item No. 22737

Overview and Properties

Synonym: Heat Shock Cognate 70

Source: N-terminal Histidine-tagged Hsc70 (human, recombinant) expressed in E. coli

Amino acids: 2-646 (full length)

Uniprot No.: P11142 Molecular Weight: 72.9 kDa

-80°C (as supplied); avoid freeze/thaw cycles and it is recommended to aliquot the Storage:

protein.

Stability: ≥1 year

≥90% estimated by SDS-PAGE **Purity:**

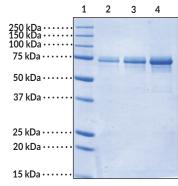
Supplied in: 50 mM HEPES, pH 8.0, and 150 mM sodium chloride

Protein

Concentration: batch specific 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: Hsc70 (1 µg) Lane 3: Hsc70 (2 µg) Lane 4: Hsc70 (4 µg)

Representative gel image shown; actual purity may vary between batches

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

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, 09/07/2021

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

Hsc70 is a chaperone from the Hsp70 chaperone family encoded by the HspA8 gene that shares 85% sequence identity to Hsp70 (encoded by HspA1A gene), and is one of the constitutively expressed members from this family. Hsc70 acts as the main housekeeping protein but is also involved in many functions such as ubiquitin-proteasome degradation pathway and transporting cytoplasmic proteins into the nucleus. Hsc70 also garnered interests as a potential therapeutic target since promoting accumulation of unfolded proteins can serve as a potential strategy to cause tumor death.

References

- 1. Stricher, F., Macri, C., and Muller, S. HSPA8/HSC70 chaperone protein: Structure, function, and chemical targeting. *Autophagy* **9(12)**, 1937-1954 (2013).
- 2. Bercovich, B., Stancovski, I., Mayer, A., et al. Ubiquitin-dependent degradation of certain protein substrates in vitro requires the molecular chaperone Hsc70. J. Biol. Chem. 272(14), 9002-9010 (1997).
- 3. Dang, C.V. and Lee, W.M. Nuclear and nucleolar targeting sequences of c-erb-A, c-myb, N-myc, p53, HSP70, and HIV tat proteins. J. Biol. Chem. 264(30), 18019-18023 (1989).
- 4. Leu, J.I., Pimkina, J., Frank, A., et al. A small molecule inhibitor of inducible heat shock protein 70. *Molecular Cell* **36(1)**, 15-27 (2009).

1180 EAST ELLSWORTH RD ANN ARBOR, MI 48108 · USA PHONE: [800] 364-9897