

# Product Information

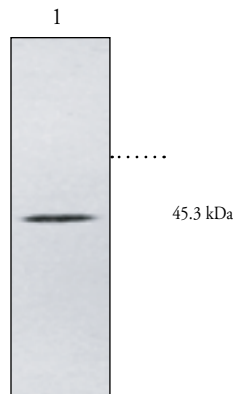


## Hepsin Polyclonal Antibody(aa 241-260)

Item No. 100022 • Lot. No. XXXXX

- Contents:** This vial contains (100-500 µg of peptide affinity-purified IgG, *lot specific*) in 500 µl TBS, pH *lot specific*, containing 50% glycerol, 0.5 mg/ml BSA, and 0.02% sodium azide
- Antigen:** Human hepsin amino acids 241-260 (GGYLPFRDPNSEENSNDIAL). The antigen alignment with other known species sequences:  
Human GGYLPFRDPNSEENSNDIAL  
Mouse GGYLPFRDP t i dNNSNDIAL  
Rat GGYLPFRDP t i dENSNDIAL
- Host:** Rabbit
- Cross-reactivity:** (+) Human hepsin; other species produced bands with unknown identity
- Storage:** ≥1 year at -20°C
- Applications:** Recommended starting dilutions: western blotting at 1 µg/ml, immunohistochemistry of formalin-fixed paraffin-embedded sections at 1-2 µg/ml. Antigen retrieval results in higher background staining and is therefore not recommended. Other applications were not attempted and therefore optimal working dilutions should be determined empirically.
- Concentration:** Varies by lot, from 0.2-1.0 mg/ml (100-500 µg/vial). Always 100 ml final working volume for western blotting.

Hepsin is a type II membrane-associated protein that has an extracellular proteolytic domain and exhibits low sequence homology to other known proteases.<sup>1,2</sup> Hepsin overexpression is observed in prostate, breast, kidney, and ovarian cancers and due to low homology to other known proteases may provide a unique target for pharmacological therapy.<sup>3-7</sup> Hepsin is necessary for cell growth *in vitro*<sup>8</sup> and may play a role in metastatic expansion by factor VII activation.<sup>6,9</sup> The observed molecular mass of hepsin by immunoblotting is 45.3 kDa. Cloning and characterization of mouse and rat hepsin indicate 88% overall homology with human hepsin.<sup>10-12</sup> Northern blotting reveals high levels of hepsin transcripts in liver.<sup>2,7</sup>



1. K562 cell lysate (25 µg)

### References

1. Tsuji, A., Torres-Rosado, A., Arai, T., *et al. J. Biol. Chem.* **266(25)**, 16948-16953 (1991).
2. Leytus, S.P., Kurachi, K., Hagen, F.S., *et al. Biochemistry* **27**, 1067-1074 (1988).
3. Dhanasekaran, S.M., Barrette, T.R., Ghosh, D., *et al. Nature* **412**, 822-826 (2001).
4. Magee, J.A., Araki, T., Patil, S., *et al. Cancer Res.* **61**, 5692-5696 (2001).
5. Welsh, J.B., Sapinoso, L.M., Su, A.I., *et al. Cancer Res.* **61**, 5974-5978 (2001).
6. Zacharski, L.R. *Thromb. Haemost.* **79**, 876-877 (1998).
7. Tanimoto, H., Yan, Y., Clarke, J., *et al. Cancer Res.* **57**, 2884-2887 (1997).
8. Torres-Rosado, A., O'Shea, K.S., Tsuji, A., *et al. Proc. Natl. Acad. Sci. USA* **90**, 7181-7185 (1993).
9. Kazama, Y., Hamamoto, T., Foster, D.C., *et al. J. Biol. Chem.* **270(1)**, 66-72 (1995).
10. Vu, T.-K.H., Liu, R.W., Haaksman, C.J., *et al. J. Biol. Chem.* **272(50)**, 31315-31320 (1997).
11. Kawamura, S., Kurachi, S., Deyashiki, Y., *et al. Eur. J. Biochem.* **262**, 755-764 (1999).
12. Farley, D., Raymond, F., and Nick, H. *Biochim. Biophys. Acta* **1173(3)**, 350-352 (1993).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/100022](http://www.caymanchem.com/catalog/100022)

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

### MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

### WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes no warranty or guarantee of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 12/15/2011

## Cayman Chemical

**Mailing address**  
1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

**Phone**  
(800) 364-9897  
(734) 971-3335

**Fax**  
(734) 971-3640

**E-Mail**  
custserv@caymanchem.com

**Web**  
[www.caymanchem.com](http://www.caymanchem.com)