

PRODUCT INFORMATION



Hormone-Sensitive Lipase (human, recombinant)

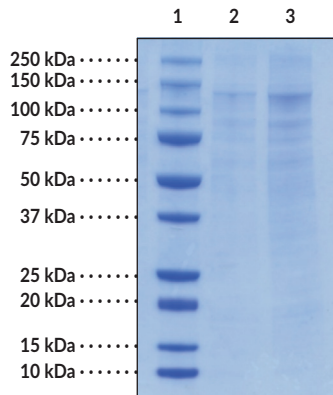
Item No. 10664

Overview and Properties

Synonyms:	HSL, LIPE
Source:	Active recombinant N-terminal hexahistidine-tagged protein, expressed in Sf21 insect cells using a baculovirus expression system.
Uniprot No.:	Q01469
Amino Acids:	2-1,076 (full-length)
Molecular Weight:	118.3 kDa
Storage:	-80°C (as supplied); avoid freeze/thaw cycles by aliquoting protein
Stability:	≥1 year
Supplied in:	50 mM sodium phosphate, 100 mM sodium chloride, 20% glycerol, pH 7.2
Protein Concentration:	<i>batch specific</i> mg/ml
Activity:	<i>batch specific</i> U/ml
Specific Activity:	<i>batch specific</i> U/mg
Unit Definition:	One unit is defined as the amount of enzyme required to produce 1 nmol of <i>p</i> -nitrophenol from <i>p</i> -nitrophenol butyrate per minute at 37°C in 100 mM sodium phosphate, pH 7.2 containing 150 mM sodium chloride and 0.5% Triton X-100

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

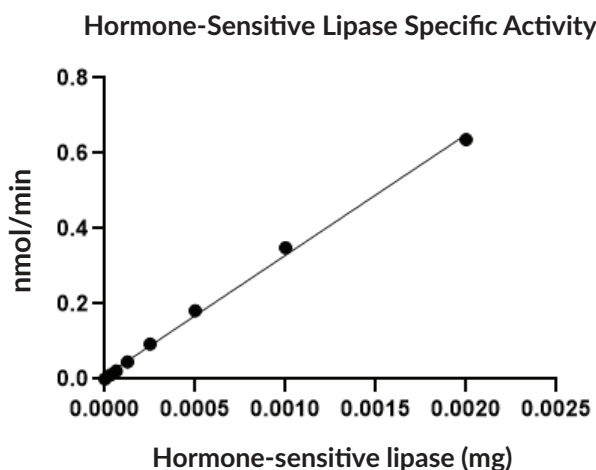
Images



Lane 1: MW Markers
Lane 2: Hormone-Sensitive Lipase (2 µg)
Lane 3: Hormone-Sensitive Lipase (4 µg)

SDS-PAGE Analysis of Hormone-Sensitive Lipase.

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, 07/17/2025

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

Hormone-sensitive lipase (HSL) is a key enzyme regulating the mobilization of fatty acids from intracellular stores.¹ There are three different isoforms of the enzyme: HSL_{adi}, which is the predominant form found in adipocytes, HSL_{tes} expressed in testes, and HSL_b found in insulin secreting β -cells.² The three isoforms are distinguished by their molecular masses.² HSL catalyses the rate limiting step in the lipolysis of triacylglycerol, found in intracellular stores, to fatty acids.³ HSL also catalyzes the hydrolysis of mono- and diacylglycerides, as well as cholesterol esters.⁴

References

1. Holm, C., Belfrage, P., Østerlund, T., et al. *Protein Eng.* **7(4)**, 537-541 (1994).
2. Krintel, C., Klint, C., Lindvall, H., et al. *PLoS One* **5(6)**, (2010).
3. Zimmermann, R., Lass, A., Haemmerle, G., et al. *Biochim. Biophys. Acta* **1791**, 494-500 (2009).
4. Saltiel, A.R. *Proc. Natl. Acad. Sci. USA* **97(2)**, 535-537 (2000).

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