

GCN5 (human), (recombinant)

hGCN5L is the catalytic histone acetyltransferase (HAT) subunit of the transcriptional coactivator complexes STAGA and TFTC.

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE272-0050	50µg
----------------	------

Manuals, SDS & CofA

[View Online »](#)

kDa

250

148

98

64

50

36

22

16



1

2

← BSA

← hGCN5

SDS-PAGE Analysis: Lane 1: MWM; Lane 2: 1.0 ug of purified GCN5 protein, with 1 mg/ml BSA.

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	GCN5L2, Histone acetyltransferase GCN5
Formulation	Liquid. In 20mM sodium citrate, pH 6.0, containing 0.25M NaCl, 1mg/ml BSA and 10% glycerol.
MW	19.4 kDa
Purity	≥95% (SDS-PAGE)
Purity Detail	Purified by multi-step chromatography.
Source	Produced in <i>E. coli</i> . Catalytic domain (aa 497-663) of human hGCN5L.
Specific Activity	≥200 pmol/min./μg assayed as production of CoA-SH from AcCoA in the presence of a peptide comprising human histone H3 residues 5-23. CoA-SH is determined with a fluorogenic, thiol-reactive maleimide (CPM).
UniProt ID	Q92830

Last modified: May 29, 2024



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)