

Lin28a (full-length) (human), (recombinant)

Lin28a is a highly conserved RNA-binding protein expressed during embryogenesis, plays roles in development, pluripotency, metabolism, tissue repair and hair regrowth.

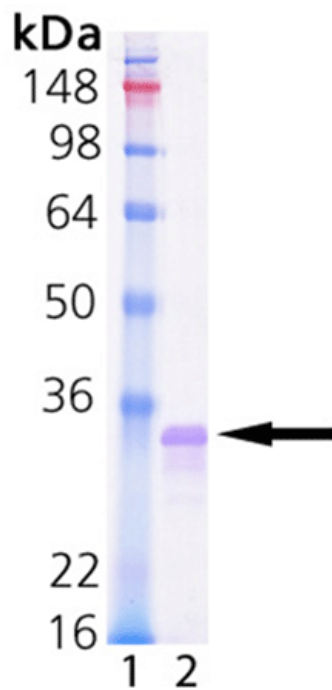
Ordering Information

[Order Online »](#)

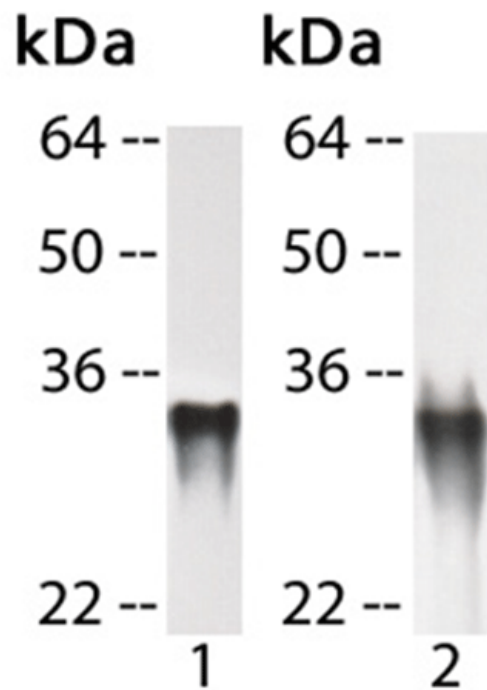
ENZ-PRT172-0050	50µg
ENZ-PRT172-0200	200µg

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis: Lane 1: MW marker; Lane 2: 1.0 μ g of purified Lin28a protein (Prod. No. ENZ-PRT172).



Western Blot analysis of 10 ng Lin28a protein (Prod. No. ENZ-PRT172) probed with Lane 1: polyclonal anti-Lin28 antibody, Lane 2: anti-6x His tag antibody.

Handling & Storage

Handling	After reconstitution, prepare aliquots and store at -80°C. Avoid freeze/thaw cycles.
Short Term Storage	-80°C
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Protein lin-28 homolog A
Formulation	Liquid. In 20mM TRIS, pH 7.2, containing 200mM sodium chloride and 5mM dithiothreitol (DTT).
MW	~24 kDa
Purity	≥95% (SDS-PAGE)
Purity Detail	Purified by multi-step chromatography.
Source	Produced in <i>E. coli</i> .
UniProt ID	Q9H9Z2

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)