

H-Ras (wild-type) (human), (recombinant)

Substrate for farnesyltransferase (>25%) with $K_m \sim 4 \mu\text{M}$ using rat brain FT.
Possesses GDP binding at >1:4 GDP/H-ras.

Citations: 2

[View Online »](#)

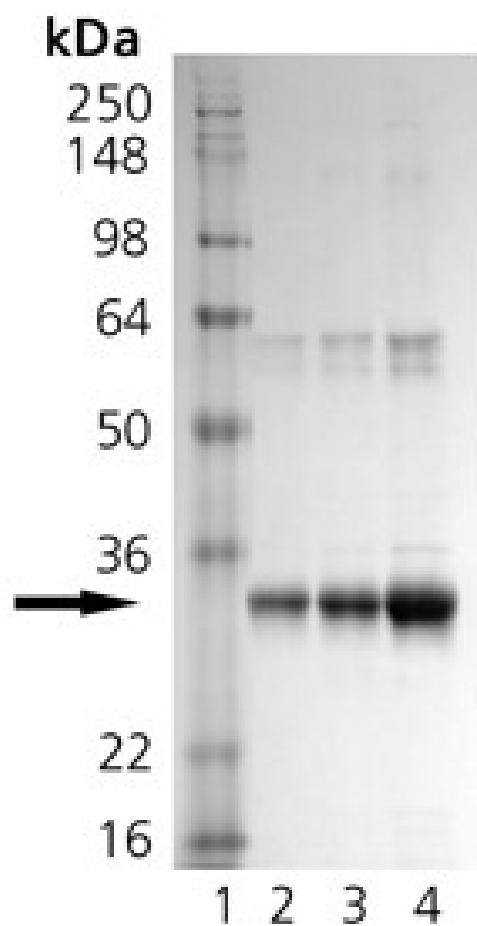
Ordering Information

[Order Online »](#)

BML-SE131-0100	100µg
----------------	-------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis: Lane 1: MW Marker, Lane 2: 1 μ g, Lane 3: 2 μ g, Lane 4: 5 μ g Ras.

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Formulation

Liquid. In 20mM TRIS-HCl, pH 8.0, containing 45mM sodium chloride, 1mM DTT, 5mM magnesium chloride and 50% glycerol.

MW

28 kDa

Purity

≥95% (SDS-PAGE)

Purity Detail

Purified by multi-step chromatography.

Source

Produced in *E. coli*.

UniProt ID

P01112

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)