

H-Ras

Harvey rat sarcoma viral oncogene homolog

Ordering Information

[Order Online »](#)

JBS-PR-107	50µg
------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Formulation liquid (Supplied in 25 mM HEPES pH 7.2, 40 mM NaCl, 3 mM DTT, 2 mM MgCl₂ and 0.01 mM GDP)

MW 21.3 kDa (189 amino acids)

Purity > 90 % (SDS-PAGE)

Source human, recombinant, E. coli

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com